

## Agilent InfinityLab LC Purification Solutions

Agilent offers tailor-made, high-efficiency LC purification solutions ranging from analytical to preparative scale. This comprehensive portfolio of state-of-the-art LC purification instruments offers you scalable LC purification solutions to meet your needs and budget. InfinityLab columns and supplies are designed to work seamlessly with Agilent InfinityLab instruments, and draw on 40+ years of technological leadership.

With Agilent OpenLab software and support from Agilent CrossLab instrument and enterprise service plans, Agilent provides a complete purification solution that will make a real difference to your lab's operational efficiency.



### **1290 Infinity II Preparative LC Systems – Purify with the new benchmark in efficiency**

Aspire to maximum efficiency in your purification workflows. Get maximum flexibility to solve all your purification challenges with highest purity, recovery, and speed.



### **1260 Infinity II Preparative LC Systems – Benefit from efficient purification – every day**

Achieve highest purity and recovery with the most affordable fraction collector. Get highest flexibility in fraction size while maintaining high-precision recovery.



### **1220/1260/1290 Infinity II Analytical-Scale LC Purification Systems – Experience the power of analytical compound isolation**

(Only preparative-scale systems are covered in this guide)

Isolate your compounds of interest through high-performance separation and low-dispersion fraction collection. Choose the bio-inert option for purification of your proteins or sensitive samples

## Pumps

|                                                                                                                                                                                 | PM*    | Part Number |
|---------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|--------|-------------|
| <b>Supplies for 1260 Infinity II Preparative Binary Pump (G7161A)</b>                                                                                                           |        |             |
| <b>Standard preventive maintenance parts</b>                                                                                                                                    |        |             |
| Preventive maintenance kit, for the 1260 Infinity II preparative binary pump (G7161A), and 1290 Infinity II preparative binary pump (G7161B) equipped with 50 mL/min pump heads | PM kit | G7161-68730 |
| Solvent bottle inlet filter, preparative, 20 µm, SST, 2/pk                                                                                                                      | 2 each | 5023-3115   |
| Filter assembly, preparative, 12.7 mm id, 2 µm, SST                                                                                                                             | 1 each | 5022-2192   |
| Piston Seal, for 1260 Infinity II preparative binary pump and 1290 Infinity II preparative binary pump G7161B with AtP pump head                                                | 4 each | 5067-4299   |
| Seal wash PE for 1260 Infinity II preparative binary pump G7161A and for AtP pump head, 1/pk                                                                                    | 4 each | 5067-6587   |
| Peristaltic pump with PharMed tubing                                                                                                                                            | 1 each | 5065-4445   |
| <b>Extended preventive maintenance parts</b>                                                                                                                                    |        |             |
| Piston collar, 5 mm, for G7161A 1260 Infinity II preparative binary pump and G7161B 1290 Infinity II preparative binary pump with AtP pump head                                 |        | G7161-60005 |
| Check valve assembly, inlet/outlet, for G7161A/B preparative pump, with 50 mL/min pump head                                                                                     |        | 5067-6642   |
| Purge valve prep                                                                                                                                                                |        | G7161-60036 |
| Solvent mixer                                                                                                                                                                   |        | G1312-87330 |
| Silicone tubing, 1 mm id, 3 mm od, 5 m                                                                                                                                          |        | 5065-9978   |
| <b>Other supplies</b>                                                                                                                                                           |        |             |
| Bottle head assembly for 1260/1290 Infinity II Preparative LC system                                                                                                            |        | G7161-60042 |
| InfinityLab capillary kit, for flow rate range 4 to 8 mL/min or column with 9.4 mm id                                                                                           |        | 5067-7015   |
| InfinityLab capillary kit, for flow rate range 15 to 40 mL/min or column with 21.2 mm id                                                                                        |        | 5067-7016   |
| InfinityLab capillary kit, for flow rate range 40 to 80 mL/min or column with 30 mm id                                                                                          |        | 5067-7017   |
| <b>InfinityLab Supplies for 1290 Infinity II Preparative Binary Pump G7161B</b>                                                                                                 |        |             |
| <b>Standard preventive maintenance parts (200 mL/min flow pump heads)</b>                                                                                                       |        |             |
| Preventive maintenance kit 1290 Infinity II Preparative Binary Pump                                                                                                             | PM kit | G7161-68740 |
| Piston seal, for 1290 Infinity II preparative binary pump G7161B with preparative pump head                                                                                     | 4 each | 5067-6589   |
| Seal wash PE for 1290 Infinity II preparative binary pump G7161B, 1/pk                                                                                                          | 4 each | 5067-6588   |
| Filter assembly, preparative, 12.7 mm id, 2 µm, SST                                                                                                                             | 1 each | 5022-2192   |
| Peristaltic pump with PharMed tubing                                                                                                                                            | 1 each | 5065-4445   |

|                                                                                                                                                                                 | PM*                     | Part Number |
|---------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|-------------------------|-------------|
| <b>Standard preventive maintenance parts (optional 50 mL pump heads)</b>                                                                                                        |                         |             |
| Preventive maintenance kit, for the 1260 Infinity II preparative binary pump (G7161A), and 1290 Infinity II preparative binary pump (G7161B) equipped with 50 mL/min pump heads | PM kit                  | G7161-68730 |
| Piston Seal, for 1260 Infinity II preparative binary pump and 1290 Infinity II preparative binary pump G7161B with AtP pump head                                                | 4 each/option depending | 5067-4299   |
| Seal wash PE for 1260 Infinity II preparative binary pump G7161A and for ATP pump head, 1/pk                                                                                    | 4 each/option depending | 5067-6587   |
| Solvent bottle inlet filter, preparative, 20 µm, SST, 2/pk                                                                                                                      | 2 each                  | 5023-3115   |
| <b>Extended preventive maintenance parts</b>                                                                                                                                    |                         |             |
| Check valve assembly double seat, inlet/outlet                                                                                                                                  |                         | G7161-60052 |
| Piston, for G7161B 1290 Infinity II preparative binary pump with preparative pump head                                                                                          |                         | G7161-60028 |
| Purge valve prep                                                                                                                                                                |                         | G7161-60036 |
| Solvent Mixer                                                                                                                                                                   |                         | G1312-87330 |
| Silicone tubing, 1 mm id, 3 mm od, 5 m                                                                                                                                          |                         | 5065-9978   |
| Piston collar, 5 mm, for G7161A 1260 Infinity II preparative binary pump and G7161B 1290 Infinity II preparative binary pump with AtP pump head                                 |                         | G7161-60005 |
| Check valve assembly, inlet/outlet, for G7161A/B preparative pump, with 50 mL/min pump head                                                                                     |                         | 5067-6642   |
| <b>Other supplies</b>                                                                                                                                                           |                         |             |
| Bottle head assembly, for 1260/1290 Infinity II Preparative LC system                                                                                                           |                         | G7161-60042 |
| InfinityLab capillary kit, for flow rate range 4 to 8 mL/min or column with 9.4 mm id                                                                                           |                         | 5067-7015   |
| InfinityLab capillary kit, for flow rate range 15 to 40 mL/min or column with 21.2 mm id                                                                                        |                         | 5067-7016   |
| InfinityLab capillary kit, for flow rate range 40 to 80 mL/min or column with 30 mm id                                                                                          |                         | 5067-7017   |
| InfinityLab capillary kit, for flow rate range 80 to 200 mL/min or column with 50 mm id                                                                                         |                         | 5067-7018   |

**\* PM (preventive maintenance) part information key**

x each: the number or amount of the part needed per PM

PM kit: a kit that includes all the standard parts for PM

optional: a part only required for PM if referring [the referenced] option is installed

## Sample injection modules

|                                                                                                                                    | PM*                     | Part Number                 |
|------------------------------------------------------------------------------------------------------------------------------------|-------------------------|-----------------------------|
| <b>InfinityLab Supplies for Preparative Open-Bed Sampler/Collector G7158B</b>                                                      |                         |                             |
| <b>Parts for preventive maintenance</b>                                                                                            |                         |                             |
| InfinityLab tubing kit, 8 mL/min, with RFID, for 1290 Infinity II preparative open-bed fraction collector                          | 1 each/option depending | <a href="#">G9321-60953</a> |
| InfinityLab tubing kit, 50 mL/min, with RFID, for 1290 Infinity II preparative open-bed fraction collector                         | 1 each/option depending | <a href="#">G9321-60952</a> |
| InfinityLab tubing kit, 150 mL/min, with RFID, for 1290 Infinity II preparative open-bed fraction collector                        | 1 each/option depending | <a href="#">G9321-60954</a> |
| InfinityLab tubing kit, 200 mL/min, with RFID, for 1290 Infinity II preparative open-bed fraction collector                        | 1 each/option depending | <a href="#">G9321-60951</a> |
| Dripping adapter, for fraction valve                                                                                               | 1 each                  | <a href="#">G9321-26600</a> |
| Rotor Seal, 2 pos/14 port preparative valve head, 600 bar                                                                          | 1 each                  | <a href="#">5068-0323</a>   |
| Peristaltic pump head, for 1290 Infinity II Preparative Open-Bed Sampler/Collector G7158B                                          | 2 each                  | <a href="#">5067-6668</a>   |
| Needle plug assembly, for use with the 1290 Infinity II Preparative Open-Bed Sampler/Collector                                     | 1 each                  | <a href="#">G9321-60036</a> |
| Tubing kit, for sampling unit                                                                                                      | 1 each                  | <a href="#">G9321-68006</a> |
| Solvent distribution tubing kit, for use with 1290 Infinity II Preparative Open-Bed Sampler/Collector.                             | 1 each                  | <a href="#">5500-1429</a>   |
| Preparative LC needle tube kit, 1.2 × 2.0 mm calibration, for use with the 1290 Infinity II Preparative Open-Bed Sampler/Collector | 1 each                  | <a href="#">G9321-68735</a> |
| Solvent bottle inlet filter, 20 µm, glass                                                                                          | 2 each                  | <a href="#">5041-2168</a>   |
| <b>Extended preventive maintenance parts</b>                                                                                       |                         |                             |
| Stator, for 2pos/14 port valve                                                                                                     | 1 each                  | <a href="#">5068-0322</a>   |
| <b>Containers and vial plates</b>                                                                                                  |                         |                             |
| InfinityLab drawer, ambient, with RFID tag                                                                                         |                         | <a href="#">G9321-60085</a> |
| InfinityLab tube container, 10 tubes, 30 × 150 mm, ambient                                                                         |                         | <a href="#">G9321-60015</a> |
| InfinityLab tube container, 18 tubes, 25 × 150 mm, ambient                                                                         |                         | <a href="#">G9321-60025</a> |
| InfinityLab tube container, 36 tubes, 16 × 150 mm, ambient                                                                         |                         | <a href="#">G9321-60129</a> |
| InfinityLab tube container, 72 tubes, 12 × 150 mm, ambient                                                                         |                         | <a href="#">G9321-60131</a> |
| InfinityLab tube container, 10 tubes, 30 × 100 mm, ambient                                                                         |                         | <a href="#">G9321-60058</a> |
| InfinityLab tube container, 18 tubes, 25 × 100 mm, ambient                                                                         |                         | <a href="#">G9321-60035</a> |
| InfinityLab tube container, 36 tubes, 16 × 100 mm, ambient                                                                         |                         | <a href="#">G9321-60055</a> |
| InfinityLab tube container, 72 tubes, 12 × 100 mm, ambient                                                                         |                         | <a href="#">G9321-60045</a> |

|                                                                                                                                          | PM* | Part Number                 |
|------------------------------------------------------------------------------------------------------------------------------------------|-----|-----------------------------|
| InfinityLab Drawer Gilson 207, ambient, contains two drawers for Gilson rack type 207, for use with 16 × 100 mm (od × length) test tubes |     | <a href="#">G9321-60201</a> |
| InfinityLab Drawer Gilson 208, ambient                                                                                                   |     | <a href="#">G9321-60208</a> |
| InfinityLab SBS well plate container, ambient                                                                                            |     | <a href="#">G9321-60051</a> |
| Vial plate, for 15 × 6 mL vials, low profile, nonstackable                                                                               |     | <a href="#">5043-1826</a>   |
| Vial plate, for 40 × 2 mL vials                                                                                                          |     | <a href="#">5023-2471</a>   |
| 96-well plate, 2.0 mL, square wells, U shape, polypropylene, 41 mm, 30/pk                                                                |     | <a href="#">5043-9300</a>   |
| Sealing mat, 96 wells, square, preslitted, silicone, 50/pk                                                                               |     | <a href="#">5043-9319</a>   |
| 96-well plate, 0.90 mL, round wells, U shape, polypropylene, 32 mm, 50/pk                                                                |     | <a href="#">5043-9305</a>   |
| 96-well plate, 1.0 mL, round wells, U shape, polypropylene, 27 mm, 50/pk                                                                 |     | <a href="#">5043-9309</a>   |
| Sealing mat, 96 wells, round, preslitted, silicone, 50/pk                                                                                |     | <a href="#">5043-9317</a>   |
| <b>Fraction tubes and vials</b>                                                                                                          |     |                             |
| Test tubes, 30 × 150 mm, 78 mL, 100/pk                                                                                                   |     | <a href="#">5190-9090</a>   |
| Test tubes, 25 × 150 mm, 55 mL, 100/pk                                                                                                   |     | <a href="#">5190-9091</a>   |
| Test tubes, 16 × 150 mm, 21 mL, 250/pk                                                                                                   |     | <a href="#">5190-9092</a>   |
| Test tubes, 12 × 150 mm, 11 mL, 250/pk                                                                                                   |     | <a href="#">5190-9093</a>   |
| Round bottom glass vial, 30 × 100 mm 100/pk                                                                                              |     | <a href="#">5042-6458</a>   |
| Round bottom glass vial, 25 × 100 mm 100/pk                                                                                              |     | <a href="#">5042-6459</a>   |
| Test tubes, 16 × 100 mm, 250/pk                                                                                                          |     | <a href="#">5022-6532</a>   |
| Test tubes, 12 × 100 mm, 250/pk                                                                                                          |     | <a href="#">5022-6531</a>   |
| Vial, screw top, clear, certified, flat bottom, 6 mL, for LC, 100/pk                                                                     |     | <a href="#">9301-1377</a>   |
| Vial, screw top, clear, high recovery, 5 mL, for LC, 30/pk                                                                               |     | <a href="#">5188-5369</a>   |
| Cap, screw, for 6 mL vials, 100/pk                                                                                                       |     | <a href="#">9301-1379</a>   |
| Septa, for 6 mL vials, 100/pk                                                                                                            |     | <a href="#">9301-1378</a>   |
| Vial, screw top, amber, write-on spot, certified, 2 mL, 100/pk                                                                           |     | <a href="#">5182-0716</a>   |
| Vial, screw top, clear, write-on spot, certified, 2 mL, 100/pk                                                                           |     | <a href="#">5182-0715</a>   |
| Cap, screw, blue, preslit PTFE/silicone, 100/pk                                                                                          |     | <a href="#">5183-2076</a>   |

|                                                                                                                                                                                 | PM*    | Part Number |
|---------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|--------|-------------|
| <b>Other supplies</b>                                                                                                                                                           |        |             |
| Sample loop, stainless steel, 400 µL, 1/16 in, for 1290 Infinity II Series Prep Open-Bed ALS/Collector (G7158B)                                                                 |        | 5068-0348   |
| Sample loop, stainless steel, 2 mL, 1/16 in                                                                                                                                     |        | 5068-0350   |
| Preparative sample loop, stainless steel, 5 mL                                                                                                                                  |        | 5068-0351   |
| Sample loop, stainless steel, 5 mL, 1/8 in, for 1290 Infinity II Series Prep Open-Bed ALS/Collector G7158B                                                                      |        | 5068-0334   |
| Sample loop, stainless steel, 10 mL, 1/8 in, for 1290 Infinity II Series Prep Open-Bed ALS/Collector                                                                            |        | 5068-0335   |
| Sample loop, stainless steel, 20 mL, 1/8 in, for 1290 Infinity II Series Prep Open-Bed ALS/Collector                                                                            |        | 5068-0336   |
| Delay and checkout calibrant, for Agilent purification systems. Two ampoules of 5 mL, Patent Blue VF sodium salt, Sunset Yellow FCF, Sudan Orange G, dissolved in DMSO, 2 mg/mL |        | 5190-8223   |
| InfinityLab bottom fume hood kit, for Agilent 1290 Infinity II preparative open-bed sampler/collector                                                                           |        | G9321-68300 |
| <b>Supplies for 1260 Infinity II Preparative Autosampler (G7157A)</b>                                                                                                           |        |             |
| <b>Parts for preventive maintenance</b>                                                                                                                                         |        |             |
| Preventive maintenance kit, for the 1260 Infinity II preparative autosampler G7157A                                                                                             | PM kit | G7157-68730 |
| Rotor seal, 6 port injection valve, preparative autosampler G7157A                                                                                                              | 1 each | 5068-0268   |
| Needle assembly, for G7157A preparative autosampler                                                                                                                             | 1 each | G7157-87201 |
| Needle seat, for G7157A preparative autosampler                                                                                                                                 | 1 each | G7157-87101 |
| Peristaltic pump with PharMed tubing                                                                                                                                            | 1 each | 5065-4445   |
| Finger caps 15/pk                                                                                                                                                               | 1 each | 5063-6506   |
| <b>Other supplies</b>                                                                                                                                                           |        |             |
| Drawer, for 66 × 2 mL vials                                                                                                                                                     | x      | G7129-60010 |
| Drawer for 18 × 6 mL vials                                                                                                                                                      | x      | G7129-60110 |
| Drawer kit for 100 × 2 mL vials, classic style                                                                                                                                  | x      | G7129-68210 |
| Multi draw loop kit, 5 mL, for G7157A, 1260 Infinity II preparative autosampler                                                                                                 | x      | G7157-68711 |
| Sample loop, stainless steel, 900 µL                                                                                                                                            | x      | G7157-60515 |
| <b>Supplies for 1260 Infinity II Manual Preparative Injector (G1328D)</b>                                                                                                       |        |             |
| <b>Parts for preventive maintenance</b>                                                                                                                                         |        |             |
| Rebuild kit, for 1260 Infinity II preparative manual injector                                                                                                                   | PM kit | 5068-0319   |
| Rotor seal for Rheodyne 3725, 3725-038, PEEK                                                                                                                                    | 1 each | 0101-1233   |
| Stator face seal, for 1260 Infinity II preparative manual injector                                                                                                              | 1 each | 5068-0321   |

|                                                                                     | PM* | Part Number |
|-------------------------------------------------------------------------------------|-----|-------------|
| <b>Other supplies</b>                                                               |     |             |
| Sample loop, stainless steel, 2 mL, 1.00 mm id                                      |     | 0101-1244   |
| Sample loop, stainless steel, 5 mL, 1.00 mm id                                      |     | 0101-1243   |
| Adapter, stainless steel, female 10 to 32 to male 5/16-24                           |     | 5067-6845   |
| Sample loop, stainless steel, 5 mL, 1/8 in od                                       |     | 5067-7031   |
| Sample loop, stainless steel, 10 mL, 1/8 in od                                      |     | 5067-7032   |
| Sample loop, stainless steel, 20 mL, 1/8 in od                                      |     | 5067-7033   |
| Sample loop, PEEK, 2 mL, 1.60 mm id                                                 |     | 0101-1229   |
| Sample loop, PEEK, 5 mL, 1.60 mm id                                                 |     | 0101-1228   |
| Sample loop, PEEK, 10 mL, 2.00 mm id                                                |     | 0101-1227   |
| Sample loop, PEEK, 20 mL, 2.00 mm id                                                |     | 0101-1226   |
| Manual syringe, 25 mL, removable Luer lock with push button valve, PTFE-tip plunger |     | 5190-1544   |
| Glass syringe, 1 mL, Luer tip, 0.05 scale                                           |     | 9301-1446   |
| Needle, sample loading, stainless steel, Luer, 16 gauge, 1/pk                       |     | RE3725086   |
| Stator manual prep injector                                                         |     | 5068-0320   |
| Holder, for mounting the preparative manual injector to the prep column organizer   |     | 5067-6831   |
| <b>Vials and Caps for use with Autosampler and Manual Injector</b>                  |     |             |
| Vial, screw top, clear, certified, flat bottom, 6 mL, for LC, 100/pk                |     | 9301-1377   |
| Vial, screw top, clear, high recovery, 5 mL, for LC, 30/pk                          |     | 5188-5369   |
| Cap, screw, for 6 mL vials, 100/pk                                                  |     | 9301-1379   |
| Septa, for 6 mL vials, 100/pk                                                       |     | 9301-1378   |
| Vial, screw top, amber, write-on spot, certified, 2 mL, 100/pk                      |     | 5182-0716   |
| Vial, screw top, clear, write-on spot, certified, 2 mL, 100/pk                      |     | 5182-0715   |
| Cap, screw, blue, preslit PTFE/silicone, 100/pk                                     |     | 5183-2076   |

**\* PM (preventive maintenance) part information key**

x each: the number or amount of the part needed per PM

PM kit: a kit that includes all the standard parts for PM

optional: a part only required for PM if referring [the referenced] option is installed

## Fraction collectors

|                                                                                                                                                                                 | PM*             | Part Number                 |
|---------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|-----------------|-----------------------------|
| <b>Supplies for 1260 Infinity II Preparative Fraction Collector (G1364E)</b>                                                                                                    |                 |                             |
| <b>Parts for preventive maintenance (G1364E)</b>                                                                                                                                |                 |                             |
| Tubing kit, for flow rate 40 to 100 mL/min, 30 mm columns, for fraction collector G1364E                                                                                        | 1 each          | <a href="#">G1364-68605</a> |
| Tubing kit, for flow rate 4 to 8 mL/min, 9.4 mm columns, for fraction collector G1364E/F and G5664B                                                                             | 1 each/optional | <a href="#">G1364-68603</a> |
| Tubing kit, for flow rate 15 to 40 mL/min, 21.2 mm columns, for fraction collector G1364E                                                                                       | 1 each/optional | <a href="#">G1364-68604</a> |
| <b>Other supplies</b>                                                                                                                                                           |                 |                             |
| Tray holding 40 tubes, 30 × 100 mm, 50 mL, for G1364E fraction collectors                                                                                                       | x               | <a href="#">G1364-84543</a> |
| Round bottom glass vial, 30 × 100 mm 100/pk                                                                                                                                     | x               | <a href="#">5042-6458</a>   |
| Tray holding 60 tubes, 25 × 100 mm, 35 mL, for G1364E fraction collectors                                                                                                       | x               | <a href="#">G1364-84544</a> |
| Round bottom glass vial, 25 × 100 mm 100/pk                                                                                                                                     | x               | <a href="#">5042-6459</a>   |
| Tray holding 126 tubes, 16 × 100 mm, 14 mL, for G1364E fraction collectors                                                                                                      | x               | <a href="#">G1364-84545</a> |
| Test tubes, 16 × 100 mm, 250/pk                                                                                                                                                 | x               | <a href="#">5022-6532</a>   |
| Tray holding 215 tubes, 12 × 100 mm, 7 mL, for G1364E fraction collectors                                                                                                       | x               | <a href="#">G1364-84536</a> |
| Test tubes, 12 × 100 mm, 250/pk                                                                                                                                                 | x               | <a href="#">5022-6531</a>   |
| Tray holding 40 × 2 mL vials, for fraction collector G1364E/F and G5664B                                                                                                        | x               | <a href="#">G1364-44612</a> |
| Tray, holding 15 × 6 mL vials, for fraction collector G1364E/F and G5664B                                                                                                       | x               | <a href="#">G1364-44613</a> |
| Tray, holding 100 × 2 mL vials, for fraction collector G1364E/F and G5664B                                                                                                      | x               | <a href="#">G1364-44610</a> |
| Tray for 100 × 2 mL vials, coolable                                                                                                                                             |                 | <a href="#">G1364-60111</a> |
| Tray, thermostatable for 4 well plates, for use with fraction collectors G1364B/F and G5664B                                                                                    | x               | <a href="#">G1364-60621</a> |
| Delay and checkout calibrant, for Agilent purification systems. Two ampoules of 5 mL, Patent Blue VF sodium salt, Sunset Yellow FCF, Sudan Orange G, dissolved in DMSO, 2 mg/mL | x               | <a href="#">5190-8223</a>   |
| Check-out solution                                                                                                                                                              | x               | <a href="#">5190-6886</a>   |
| Back pressure regulator, 20 psi                                                                                                                                                 | x               | <a href="#">5067-6840</a>   |
| Exhaust tube adapter kit, for Agilent 1260 Infinity II Preparative/Analytical fraction collector, use with G1364E/F, G5664B                                                     |                 | <a href="#">G1364-68709</a> |

|                                                                                                                                          | PM*                     | Part Number                 |
|------------------------------------------------------------------------------------------------------------------------------------------|-------------------------|-----------------------------|
| <b>InfinityLab Supplies for 1290 Infinity II Preparative OpenBed Fraction Collector (G7159B)</b>                                         |                         |                             |
| <b>Parts for preventive maintenance</b>                                                                                                  |                         |                             |
| InfinityLab tubing kit, 8 mL/min, with RFID, for 1290 Infinity II preparative open-bed fraction collector                                | 1 each/option depending | <a href="#">G9321-60953</a> |
| InfinityLab tubing kit, 50 mL/min, with RFID, for 1290 Infinity II preparative open-bed fraction collector                               | 1 each/option depending | <a href="#">G9321-60952</a> |
| InfinityLab tubing kit, 150 mL/min, with RFID, for 1290 Infinity II preparative open-bed fraction collector                              | 1 each/option depending | <a href="#">G9321-60954</a> |
| InfinityLab tubing kit, 200 mL/min, with RFID, for 1290 Infinity II preparative open-bed fraction collector                              | 1 each/option depending | <a href="#">G9321-60951</a> |
| Dripping adapter, for fraction valve                                                                                                     | 1 each                  | <a href="#">G9321-26600</a> |
| <b>Containers and vial plates</b>                                                                                                        |                         |                             |
| InfinityLab drawer, ambient, with RFID tag                                                                                               |                         | <a href="#">G9321-60085</a> |
| InfinityLab tube container, 10 tubes, 30 × 150 mm, ambient                                                                               |                         | <a href="#">G9321-60015</a> |
| InfinityLab tube container, 18 tubes, 25 × 150 mm, ambient                                                                               |                         | <a href="#">G9321-60025</a> |
| InfinityLab tube container, 36 tubes, 16 × 150 mm, ambient                                                                               |                         | <a href="#">G9321-60129</a> |
| InfinityLab tube container, 72 tubes, 12 × 150 mm, ambient                                                                               |                         | <a href="#">G9321-60131</a> |
| InfinityLab tube container, 10 tubes, 30 × 100 mm, ambient                                                                               |                         | <a href="#">G9321-60058</a> |
| InfinityLab tube container, 18 tubes, 25 × 100 mm, ambient                                                                               |                         | <a href="#">G9321-60035</a> |
| InfinityLab tube container, 36 tubes, 16 × 100 mm, ambient                                                                               |                         | <a href="#">G9321-60055</a> |
| InfinityLab tube container, 72 tubes, 12 × 100 mm, ambient                                                                               |                         | <a href="#">G9321-60045</a> |
| InfinityLab Drawer Gilson 207, ambient, contains two drawers for Gilson rack type 207, for use with 16 × 100 mm (od × length) test tubes |                         | <a href="#">G9321-60201</a> |
| InfinityLab Drawer Gilson 208, ambient                                                                                                   |                         | <a href="#">G9321-60208</a> |
| InfinityLab SBS well plate container, ambient                                                                                            |                         | <a href="#">G9321-60051</a> |
| Vial plate, for 15 × 6 mL vials, low profile, nonstackable                                                                               |                         | <a href="#">5043-1826</a>   |
| Vial plate, for 40 × 2 mL vials                                                                                                          |                         | <a href="#">5023-2471</a>   |
| 96-well plate, 2.0 mL, square wells, U shape, polypropylene, 41 mm, 30/pk                                                                |                         | <a href="#">5043-9300</a>   |
| Sealing mat, 96 wells, square, preslitted, silicone, 50/pk                                                                               |                         | <a href="#">5043-9319</a>   |
| 96-well plate, 0.90 mL, round wells, U shape, polypropylene, 32 mm, 50/pk                                                                |                         | <a href="#">5043-9305</a>   |

|                                                                                                                                                                                 | PM* | Part Number |
|---------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|-----|-------------|
| 96-well plate, 1.0 mL, round wells, U shape, polypropylene, 27 mm, 50/pk                                                                                                        |     | 5043-9309   |
| Sealing mat, 96 wells, round, preslitted, silicone, 50/pk                                                                                                                       |     | 5043-9317   |
| <b>Fraction tubes and vials</b>                                                                                                                                                 |     |             |
| Test tubes, 30 × 150 mm, 78 mL, 100/pk                                                                                                                                          |     | 5190-9090   |
| Test tubes, 25 × 150 mm, 55 mL, 100/pk                                                                                                                                          |     | 5190-9091   |
| Test tubes, 16 × 150 mm, 21 mL, 250/pk                                                                                                                                          |     | 5190-9092   |
| Test tubes, 12 × 150 mm, 11 mL, 250/pk                                                                                                                                          |     | 5190-9093   |
| Round bottom glass vial, 30 × 100 mm 100/pk                                                                                                                                     |     | 5042-6458   |
| Round bottom glass vial, 25 × 100 mm 100/pk                                                                                                                                     |     | 5042-6459   |
| Test tubes, 16 × 100 mm, 250/pk                                                                                                                                                 |     | 5022-6532   |
| Test tubes, 12 × 100 mm, 250/pk                                                                                                                                                 |     | 5022-6531   |
| Vial, screw top, clear, certified, flat bottom, 6 mL, for LC, 100/pk                                                                                                            |     | 9301-1377   |
| Vial, screw top, clear, high recovery, 5 mL, for LC, 30/pk                                                                                                                      |     | 5188-5369   |
| Cap, screw, for 6 mL vials, 100/pk                                                                                                                                              |     | 9301-1379   |
| Septa, for 6 mL vials, 100/pk                                                                                                                                                   |     | 9301-1378   |
| Vial, screw top, amber, write-on spot, certified, 2 mL, 100/pk                                                                                                                  |     | 5182-0716   |
| Vial, screw top, clear, write-on spot, certified, 2 mL, 100/pk                                                                                                                  |     | 5182-0715   |
| Cap, screw, blue, preslit PTFE/silicone, 100/pk                                                                                                                                 |     | 5183-2076   |
| <b>Other supplies</b>                                                                                                                                                           |     |             |
| 20 psi Backpressure regulator                                                                                                                                                   |     | 5067-6840   |
| Delay and checkout calibrant, for Agilent purification systems. Two ampoules of 5 mL, Patent Blue VF sodium salt, Sunset Yellow FCF, Sudan Orange G, dissolved in DMSO, 2 mg/mL |     | 5190-8223   |
| <b>Supplies for 1260 Infinity II Preparative Valve-Based Fraction Collector (G7166A)</b>                                                                                        |     |             |
| Delay and checkout calibrant, for Agilent purification systems. Two ampoules of 5 mL, Patent Blue VF sodium salt, Sunset Yellow FCF, Sudan Orange G, dissolved in DMSO, 2 mg/mL |     | 5190-8223   |
| InfinityLab solvent bottle, clear 125 mL, with cap                                                                                                                              |     | 9301-6527   |
| InfinityLab solvent bottle, clear, 500 mL, with cap                                                                                                                             |     | 9301-6523   |
| InfinityLab solvent bottle, clear, 1,000 mL, with cap                                                                                                                           |     | 9301-6528   |
| InfinityLab identification silicone rings, 4 colors, 8/pk                                                                                                                       |     | 9301-6529   |
| InfinityLab stickers, for solvent bottles, 100/pk                                                                                                                               |     | 9301-6530   |
| InfinityLab fitting, for 2.3 mm od tubing, 2/pk                                                                                                                                 |     | 5043-1215   |

## UV detectors

|                                                                                                                                               | PM*    | Part Number |
|-----------------------------------------------------------------------------------------------------------------------------------------------|--------|-------------|
| <b>1260 Infinity II Diode Array &amp; Multi Wavelength Detector (G7115A/G7165A)</b>                                                           |        |             |
| InfinityLab long-life deuterium DAD/MWD lamp, with RFID tag                                                                                   | 1 each | 2140-0820   |
| Tungsten VIS lamp                                                                                                                             |        | G1103-60001 |
| InfinityLab preparative flow cell, for MWD (G7165A) and DAD (G7115A), 3 mm path length                                                        |        | G7115-60001 |
| InfinityLab preparative flow cell, for MWD (G7165A) and DAD (G7115A), 0.3 mm path length                                                      |        | G7115-60002 |
| InfinityLab preparative flow cell, for MWD (G7165A) and DAD (G7115A) 0.06 mm path length                                                      |        | G7115-60003 |
| Capillary, stainless steel, for preparative flow cells (DAD/MWD), use for flow rates 4 to 8 mL/min. Two capillaries required per flow cell.   |        | 5500-1342   |
| Capillary, stainless steel, for preparative flow cells (DAD/MWD), use for flow rates 15 to 40 mL/min. Two capillaries required per flow cell. |        | 5500-1343   |
| Tubing, PTFE-ESD, 80 cm length, 0.5 mm id, 1.6 mm od, with fittings for preparative flow cell                                                 |        | G1315-67306 |
| <b>1260 Infinity II Variable Wavelength Detector (G7114A)</b>                                                                                 |        |             |
| InfinityLab long-life deuterium VWD lamp, with RFID tag                                                                                       | 1 each | G1314-60101 |
| Preparative flow cell for VWD, 3 mm path length, 2 µl cell volume, with RFID tag                                                              |        | G1314-60024 |
| Preparative flow cell for VWD, 0.3 mm path length, with RFID tag                                                                              |        | G1314-60025 |
| Preparative flow cell for VWD, 0.06 mm path length, with RFID tag                                                                             |        | G1314-60023 |
| Preparative flow cell, Quartz, 3 mm, for VWD                                                                                                  |        | G7114-60024 |
| <b>Supplies for 1260 Infinity II Evaporative Light Scattering Detector</b>                                                                    |        |             |
| Splitter kit, 2 to 10 mL/min, analytical, for Agilent G4260B evaporative light scattering detector                                            |        | 5023-2256   |
| Splitter kit, 8 to 40 mL/min, preparative, for Agilent G4260B evaporative light scattering detector                                           |        | 5023-2257   |

### \* PM (preventive maintenance) part information key

x each: the number or amount of the part needed per PM

PM kit: a kit that includes all the standard parts for PM

optional: a part only required for PM if referring [the referenced] option is installed

## Column organizers

|                                                                                        | PM* | Part Number                 |
|----------------------------------------------------------------------------------------|-----|-----------------------------|
| <b>Supplies for Preparative Column Compartment (G7163B) /Column Organizer (G9328A)</b> |     |                             |
| Column holder clips (lamella type), 2/pk                                               |     | <a href="#">G7116-68003</a> |
| Column holder clamp for columns 9.4 to 30 mm id                                        |     | <a href="#">5067-6849</a>   |
| Column holder clamp, for columns 30 to 50 mm id                                        |     | <a href="#">G7163-60002</a> |

## Supporting modules

|                                                                                                                               | PM*    | Part Number               |
|-------------------------------------------------------------------------------------------------------------------------------|--------|---------------------------|
| <b>Supplies for 1290 Infinity II MS Flow Modulator (G7170B)</b>                                                               |        |                           |
| Rebuild Kit, for flow modulator                                                                                               | 1 each | <a href="#">1535-5215</a> |
| InfinityLab capillary kit for mass-based fraction collection systems                                                          |        | <a href="#">5067-7023</a> |
| <b>Supplies for 1260 Infinity II Delay Coil Organizer (G9324A)</b>                                                            |        |                           |
| InfinityLab knitted delay coils kit, for use in mass-based prep systems, for flow rate 4 to 8 mL/min or 9.4 mm id columns     |        | <a href="#">5067-6800</a> |
| InfinityLab knitted delay coils kit, for use in mass-based prep systems, for flow rates 15 to 40 mL/min or 21.2 mm id columns |        | <a href="#">5067-6801</a> |
| InfinityLab knitted delay coils kit, for use in mass-based preparative systems, for flow rate 40 to 80 mL/min                 |        | <a href="#">5067-6802</a> |
| InfinityLab knitted delay coils kit, for use in mass-based preparative systems                                                |        | <a href="#">5067-6804</a> |

\* PM (preventive maintenance) part information key

x each: the number or amount of the part needed per PM

PM kit: a kit that includes all the standard parts for PM

optional: a part only required for PM if referring [the referenced] option is installed

## Valve supplies

|                                                                                                 | PM*    | Part Number               |
|-------------------------------------------------------------------------------------------------|--------|---------------------------|
| <b>Supplies for 2Pos/10Port Valve Head Prep Scale (G4730A)</b>                                  |        |                           |
| Rotor seal, five-groove, 600 bar, for 2pos/10 port preparative scale valve head (p/n 5067-4193) | 1 each | <a href="#">5068-0153</a> |
| <b>Supplies for 8Pos/9Port Valve Head Prep Scale (G4731A)</b>                                   |        |                           |
| Rotor seal, single-groove, 600 bar, 2 pos/10 port preparative scale valve head (p/n 5067-4193)  | 1 each | <a href="#">5068-0155</a> |
| <b>Supplies for 6-Column Selection Valve Head Prep Scale (G4734A), obsolete</b>                 |        |                           |
| Rotor seal, for the 6-column selector valve head, Preparative scale, 600 bar                    | 1 each | <a href="#">5068-0249</a> |

|                                                                                                                                                             | PM*    | Part Number                 |
|-------------------------------------------------------------------------------------------------------------------------------------------------------------|--------|-----------------------------|
| <b>Supplies for 6-Column Selection Valve Head Prep Scale (G4734B)</b>                                                                                       |        |                             |
| Rotor, for use with 6-column selector valve head, preparative scale (p/n <a href="#">5067-6722</a> )                                                        | 1 each | <a href="#">5068-0331</a>   |
| Bearing ring, for 2D-LC valve head, 1,300 bar                                                                                                               |        | <a href="#">1535-4045</a>   |
| Capillary kit, for Preparative column selection valve, flow rate 4 to 8 mL/min                                                                              |        | <a href="#">5067-6867</a>   |
| Capillary kit, for Preparative column selection valve, flow rate 15 to 40 mL/min                                                                            |        | <a href="#">5067-6868</a>   |
| Capillary kit, for Preparative column selection valve, flow rates 40 to 80 mL/min                                                                           |        | <a href="#">5067-6869</a>   |
| Capillary kit, for Preparative column selection valve, flow rates 80 to 200 mL/min                                                                          |        | <a href="#">5067-6870</a>   |
| <b>Supplies for 1260 Infinity II Clustering Valve (G9322A)</b>                                                                                              |        |                             |
| Tubing, PTFE (ESD), 0.8 mm id, 1.6 mm od, 6 m                                                                                                               |        | <a href="#">5043-1540</a>   |
| Capillary kit, for clustering up to 3 fraction collectors                                                                                                   |        | <a href="#">5067-6811</a>   |
| Rail system, for fixation of clustering valve G9322A or recovery valve G7166A, on Agilent 1290 Infinity II preparative open-bed fraction collector G7159A/B |        | <a href="#">G9321-60013</a> |
| Tubing guide assembly, for Agilent 1290 Infinity II preparative open-bed fraction collector G7159A/B                                                        |        | <a href="#">G9321-60026</a> |
| <b>Supplies for 2Pos/14Port Valve Head Prep (G4738A)</b>                                                                                                    |        |                             |
| Rotor Seal, 2 pos/14 port preparative valve head, 600 bar                                                                                                   | 1 each | <a href="#">5068-0323</a>   |
| InfinityLab upgrade capillary kit, for Autoscale Fraction Collection                                                                                        |        | <a href="#">5067-7019</a>   |
| <b>Miscellaneous InfinityLab supplies</b>                                                                                                                   |        |                             |
| System tool kit                                                                                                                                             |        | <a href="#">G7120-68708</a> |
| Preparative pressure relief valve kit                                                                                                                       |        | <a href="#">5067-6857</a>   |
| Back Pressure Regulator, 100 psi                                                                                                                            |        | <a href="#">5042-6443</a>   |
| Back pressure regulator, 20 psi                                                                                                                             |        | <a href="#">5067-6840</a>   |
| <b>Stay Safe Caps and Waste Management</b>                                                                                                                  |        |                             |
| InfinityLab Stay Safe cap GL45, 1 port for preparative LC                                                                                                   |        | <a href="#">5043-1333</a>   |
| InfinityLab Stay Safe cap GL45, 2 ports for preparative LC                                                                                                  |        | <a href="#">5043-1334</a>   |
| InfinityLab thread adapter, GL45(M) to GPI 38(F), PTFE                                                                                                      |        | <a href="#">5043-1192</a>   |
| Fitting for 4.7 mm tubing, for prep LC Stay Safe caps, 2/pk                                                                                                 |        | <a href="#">5043-1335</a>   |
| 2 Screw plugs for Stay Safe caps prep LC                                                                                                                    |        | <a href="#">5005-2970</a>   |
| Kit, including InfinityLab Stay Safe cap (S60 thread, 4 ports) and 10 L waste can                                                                           |        | <a href="#">5043-1338</a>   |
| Waste can, 10 L, with S60 thread                                                                                                                            |        | <a href="#">5043-1337</a>   |

|                                                                                                 | PM* | Part Number                 |
|-------------------------------------------------------------------------------------------------|-----|-----------------------------|
| <b>Tubing and Adapters</b>                                                                      |     |                             |
| Tubing, PTFE (ESD), 1.6 mm od, 0.8 mm id, 6 m                                                   |     | <a href="#">5043-1540</a>   |
| Tubing, PTFE (ESD), 1.6 mm od, 0.5 mm id, 80 cm length, with fittings for preparative flow cell |     | <a href="#">G1315-67306</a> |
| Tubing, PTFE, ESD-striped, 2.0 mm od, 1.2 mm id, 2 m length                                     |     | <a href="#">5023-2878</a>   |
| Tubing, PTFE, ESD-striped, 2.5 mm od, 1.6 mm id, 6 m length                                     |     | <a href="#">5023-2882</a>   |
| Fitting, 1/4-28, for ESD-PEEK tubing with 1.6 mm OD, 6/pk                                       |     | <a href="#">5023-2898</a>   |
| Fitting, 1/4-28, for ESD-PEEK tubing with 2.0 mm OD, 6/pk                                       |     | <a href="#">5023-2899</a>   |
| Fitting, 1/4-28, for ESD-PEEK tubing with 2.5 mm OD, 6/pk                                       |     | <a href="#">5023-2883</a>   |
| Fitting, for ESD-PEEK tubing, 5/16 to 3/16 inch, 12/pk                                          |     | <a href="#">5043-1847</a>   |
| Union, flat-bottom, 1/4-28 ESD                                                                  |     | <a href="#">5023-2508</a>   |
| Fitting adapter, 1-4 28 to 10-32 coned, ESD protective material                                 |     | <a href="#">5500-1410</a>   |
| Adapter nut, flat bottom, 1/4-28-M, 5/16-24-F, stainless steel                                  |     | <a href="#">5023-3093</a>   |
| Adapter, stainless steel, female 10-32 to male 5/16-24                                          |     | <a href="#">5067-6845</a>   |
| Adapter, stainless steel, female to male 10-32 to 1/4-28                                        |     | <a href="#">5023-1803</a>   |
| Adapter, stainless steel, 10-32 female to 1/4-28 female, for prep LC                            |     | <a href="#">5023-2875</a>   |
| Adapter, stainless steel, female 1/4-28 to male 10-32, for prep LC                              |     | <a href="#">5023-2512</a>   |
| <b>System Capillaries</b>                                                                       |     |                             |
| Capillary, stainless steel, 0.12 × 70 mm, SI/SI                                                 |     | <a href="#">5500-1329</a>   |
| Capillary, stainless steel, 0.12 × 300 mm, NX/NX                                                |     | <a href="#">5500-1432</a>   |
| Capillary, stainless steel, 12 × 350 mm, SX/U6                                                  |     | <a href="#">5500-1433</a>   |
| Capillary, stainless steel, 0.12 × 2000 mm, SX/U6                                               |     | <a href="#">5500-1434</a>   |
| Capillary, stainless steel, 0.17 × 800 mm, NX/NX                                                |     | <a href="#">5500-1431</a>   |
| Capillary, stainless steel, 0.17 × 800 mm, SX                                                   |     | <a href="#">5500-1430</a>   |
| Capillary, stainless steel, 0.17 × 1150 mm, SX                                                  |     | <a href="#">5500-1423</a>   |
| Capillary, stainless steel, 0.17 × 1200 mm, NX/NX                                               |     | <a href="#">5500-1462</a>   |
| Capillary, stainless steel, 0.25 × 220 mm, NX/NX                                                |     | <a href="#">5500-1425</a>   |
| Capillary, stainless steel, 0.25 × 700 mm, SI/SI                                                |     | <a href="#">5500-1330</a>   |
| Capillary, stainless steel, 0.25 × 800 mm, NX/NX                                                |     | <a href="#">5500-1461</a>   |
| Capillary, stainless steel, 0.25 × 1100 mm, NX/NX                                               |     | <a href="#">5500-1422</a>   |
| Capillary, stainless steel, 0.25 × 1200 mm, NX/NX                                               |     | <a href="#">5500-1424</a>   |
| Capillary, stainless steel, 1.0 × 1900 mm, SX/SI                                                |     | <a href="#">5500-1415</a>   |

|                                                                    | PM* | Part Number               |
|--------------------------------------------------------------------|-----|---------------------------|
| <b>0.3 mm id capillaries</b>                                       |     |                           |
| Capillary, stainless steel, 0.3 × 500 mm, SX/-                     |     | <a href="#">5500-1316</a> |
| Capillary, stainless steel, 0.3 × 500 mm, long socket, no fittings |     | <a href="#">5500-1448</a> |
| Capillary, stainless steel, 0.3 × 550 mm, SX/-                     |     | <a href="#">5500-1449</a> |
| Capillary, stainless steel, 0.3 × 550 mm, SX/-/SI                  |     | <a href="#">5500-1408</a> |
| Capillary, stainless steel, 0.3 × 650 mm, SX/-                     |     | <a href="#">5500-1314</a> |
| Capillary, stainless steel, 0.3 × 650 mm, SX/-/SX                  |     | <a href="#">5500-1308</a> |
| Capillary, stainless steel, 0.3 × 750 mm, SX/SI                    |     | <a href="#">5500-1437</a> |
| Capillary, stainless steel, 0.3 × 800 mm, SX/-                     |     | <a href="#">5500-1406</a> |
| Capillary, stainless steel, 0.3 × 900 mm, no fittings              |     | <a href="#">5500-1450</a> |
| Capillary, stainless steel, 0.3 × 1000 mm, SX/-                    |     | <a href="#">5500-1407</a> |
| Capillary, stainless steel, 0.3 × 1150 mm, SX                      |     | <a href="#">5500-1435</a> |
| Capillary, stainless steel, 0.3 × 1450 mm, NX/NX                   |     | <a href="#">5500-1436</a> |
| Capillary, stainless steel, 0.3 × 1900 mm, SX/SI                   |     | <a href="#">5500-1412</a> |
| <b>0.5 mm id capillaries</b>                                       |     |                           |
| Capillary, stainless steel, 0.5 × 150 mm, no fittings              |     | <a href="#">5022-6509</a> |
| Capillary, stainless steel, 0.5 × 280 mm, no fittings              |     | <a href="#">5500-1336</a> |
| Capillary, stainless steel, 0.5 × 400 mm, no fittings              |     | <a href="#">5500-1334</a> |
| Capillary, stainless steel, 0.5 × 500 mm, no fittings              |     | <a href="#">5500-1451</a> |
| Capillary, stainless steel, 0.5 × 500 mm, SX/-                     |     | <a href="#">5500-1317</a> |
| Capillary, stainless steel, 0.5 × 550 mm, SX/-/SI                  |     | <a href="#">5500-1386</a> |
| Capillary, stainless steel, 0.5 × 550 mm, SX/-                     |     | <a href="#">5500-1452</a> |
| Capillary, stainless steel, 0.5 × 600 mm, no fittings              |     | <a href="#">5500-1333</a> |
| Capillary, stainless steel, 0.5 × 650 mm, SX/-                     |     | <a href="#">5500-1315</a> |
| Capillary, stainless steel, 0.5 × 650 mm, SX/-/SX                  |     | <a href="#">5500-1309</a> |
| Capillary, stainless steel, 0.5 × 750 mm, SX/SI                    |     | <a href="#">5500-1440</a> |
| Capillary, stainless steel, 0.5 × 800 mm, no fittings              |     | <a href="#">5500-1335</a> |
| Capillary, stainless steel, 0.5 × 800 mm, SX/-                     |     | <a href="#">5500-1384</a> |
| Capillary, stainless steel, 0.5 × 900 mm, no fittings              |     | <a href="#">5500-1453</a> |
| Capillary, stainless steel, 0.5 × 1000 mm, SX/-                    |     | <a href="#">5500-1385</a> |
| Capillary, stainless steel, 0.5 × 1150 mm, SX/-                    |     | <a href="#">5500-1438</a> |
| Capillary, stainless steel, 0.5 × 1450 mm, NX/NX                   |     | <a href="#">5500-1439</a> |
| <b>0.6 mm id capillaries</b>                                       |     |                           |
| Capillary, stainless steel, 0.6 × 100 mm, SX/SX                    |     | <a href="#">5500-1427</a> |
| Capillary, stainless steel, 0.6 × 500 mm, SX/-                     |     | <a href="#">5500-1320</a> |
| Capillary, stainless steel, 0.6 × 500 mm, no fittings              |     | <a href="#">5500-1454</a> |
| Capillary, stainless steel, 0.6 × 550 mm, SX/-                     |     | <a href="#">5500-1455</a> |
| Capillary, stainless steel, 0.6 × 650 mm, SX/-/SX                  |     | <a href="#">5500-1318</a> |
| Capillary, stainless steel, 0.6 × 650 mm, SX/-                     |     | <a href="#">5500-1319</a> |
| Capillary, stainless steel, 0.6 × 800 mm, SX/-                     |     | <a href="#">5500-1387</a> |
| Capillary, stainless steel, 0.6 × 900 mm, no fittings              |     | <a href="#">5500-1456</a> |
| Capillary, stainless steel, 0.6 × 1000 mm, SX/-                    |     | <a href="#">5500-1388</a> |
| Capillary, stainless steel, 0.6 × 1450 mm, NX/NX                   |     | <a href="#">5500-1442</a> |
| Capillary, stainless steel, 0.6 × 1900 mm, SX/SI                   |     | <a href="#">5500-1413</a> |

|                                                           | PM* | Part Number               |
|-----------------------------------------------------------|-----|---------------------------|
| <b>0.7 mm id capillaries</b>                              |     |                           |
| Capillary, stainless steel, 0.7 × 550 mm, SX/-/SI         |     | <a href="#">5500-1389</a> |
| Capillary, stainless steel, 0.7 × 750 mm, SX/SI           |     | <a href="#">5500-1443</a> |
| Capillary, stainless steel, 0.7 × 1900 mm, SX/SI          |     | <a href="#">5500-1414</a> |
| <b>0.94 mm id capillaries</b>                             |     |                           |
| Capillary, stainless steel, 0.94 × 60 mm, SI/-            |     | <a href="#">5500-1447</a> |
| Capillary, stainless steel, 0.94 × 500 mm,<br>no fittings |     | <a href="#">5500-1457</a> |
| Capillary, stainless steel, 0.94 × 500 mm, SX/-           |     | <a href="#">5500-1391</a> |
| Capillary, stainless steel, 0.94 × 550 mm, SX/-           |     | <a href="#">5500-1458</a> |
| Capillary, stainless steel, 0.94 × 550 mm, SX/-/SI        |     | <a href="#">5500-1395</a> |
| Capillary, stainless steel, 0.94 × 650 mm, SX/-           |     | <a href="#">5500-1393</a> |
| Capillary, stainless steel, 0.94 × 650 mm, SX/-/SX        |     | <a href="#">5500-1390</a> |
| Capillary, stainless steel, 0.94 × 750 mm, SX/SI          |     | <a href="#">5500-1446</a> |
| Capillary, stainless steel, 0.94 × 800 mm, SX/-           |     | <a href="#">5500-1392</a> |
| Capillary, stainless steel, 0.94 × 900 mm,<br>no fittings |     | <a href="#">5500-1459</a> |
| Capillary, stainless steel, 0.94 × 1000 mm, SX/-          |     | <a href="#">5500-1394</a> |
| Capillary, stainless steel, 0.94 × 1150 mm, SX/-          |     | <a href="#">5500-1444</a> |
| Capillary, stainless steel, 0.94 × 1450 mm, NX/NX         |     | <a href="#">5500-1445</a> |

#### Capillary fitting codes key

SI fitting = Stainless steel fitting with short screw

SX fitting = Stainless steel fitting with extra long screw

U6 = Stainless steel female 10-32 port thread size connector

NX = Stainless steel fitting with long screw with SW5 hex head

[www.agilent.com/chem/infinitylab-lc-purification](http://www.agilent.com/chem/infinitylab-lc-purification)

DE-016397

This information is subject to change without notice.

© Agilent Technologies, Inc. 2020-2026  
Printed in the USA, July 21, 2026  
5994-2810EN