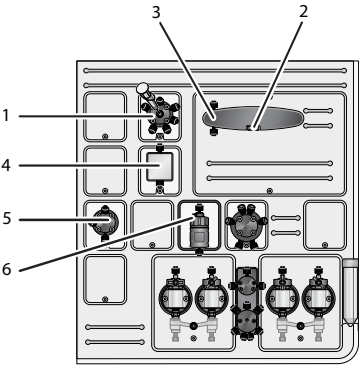


Micro kit Quick Guide

For detailed information on how to install the Micro kit refer to the *Micro kit Instructions for Use* (29337720). The quick guide provides you with an overview of the tubing used in the Blue tubing kit and the Red tubing kit.

The illustration below shows the placement of micro kit hardware modules.

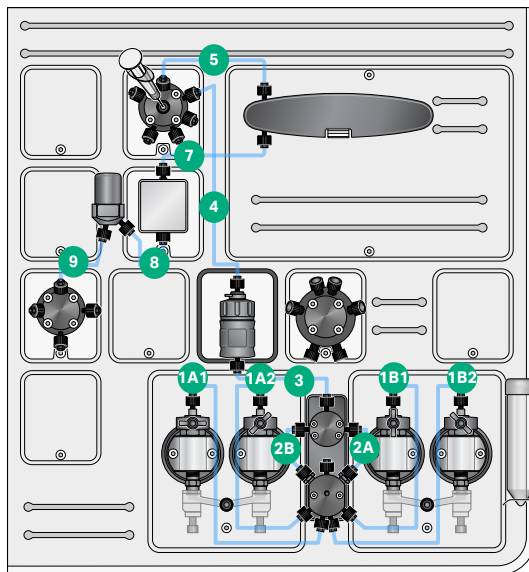


Label	Description
1	Injection valve V9M-J
2	UV monitor U9-M
3	UV flow cell U9-2M
4	Conductivity monitor C9M
5	Outlet valve V9M-Os
6	Mixer M9 with 0.6 mL or on-line filter

The table below states the information of the delay volumes with standard configuration of the Micro kit.

Tubing kit	Delay volume, Monitor to outlet valve	Delay volume, Monitor to frac F9-R or F9-T
Blue i.d. 0.25 mm	38 μ L	60 μ L
Red i.d. 0.13 mm	12 μ L	18 μ L

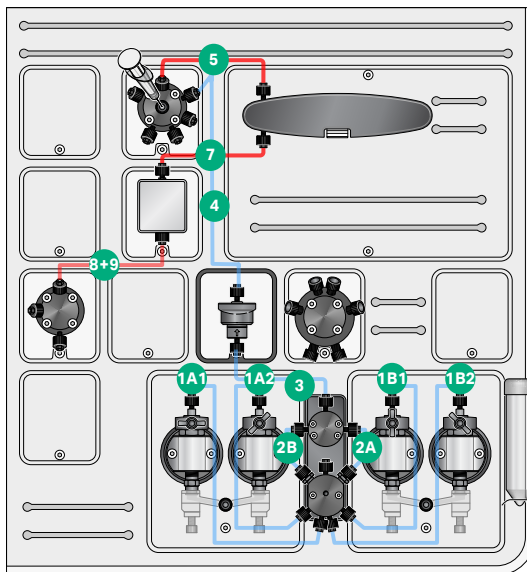
The illustration below shows the Blue tubing kit.



The table below states the information of the Blue tubing kit i.d. 0.25 mm (post-injection) for flow rates 0.25 – 2 ml/min.

Label	Tubing	Length	Description
1A1	PEEK I.D. 0.25 mm	340 mm	Pump head to pump restrictor (left side)
1A2	PEEK I.D. 0.25 mm	340 mm	Pump head to pump restrictor (left side)
1B1	PEEK I.D. 0.25 mm	340 mm	Pump head to pump restrictor (right side)
1B2	PEEK I.D. 0.25 mm	340 mm	Pump head to pump restrictor (right side)
2A	PEEK I.D. 0.25 mm	115 mm	Pump restrictor (right side) to Pressure monitor (right side)
2B	PEEK I.D. 0.25 mm	115 mm	Pump restrictor (right side) to Pressure monitor (left side)
3	PEEK I.D. 0.25 mm	160 mm	Pressure monitor to Mixer 0.6mL
4	PEEK I.D. 0.25 mm	250 mm	Mixer 0.6 mL to Injection valve V9M-J (port Pump)
5	PEEK I.D. 0.25 mm	520 mm	Injection valve V9M-J (port Col) to Column
(6)	M-M connector	N/A	Column to UV monitor U9-M
7	PEEK I.D. 0.25 mm	170 mm	UV monitor U9-M to Conductivity monitor C9M
8	PEEK I.D. 0.25 mm	95 mm	Conductivity monitor C9M to Flow restrictor (port IN)
9	PEEK I.D. 0.25 mm	135 mm	Flow restrictor to Outlet valve V9M-Os (port In)
Frac	PEEK I.D. 0.25 mm	400 mm	Outlet valve V9M-Os (port Frac) to Fraction collector nozzle

The illustration below shows the Red tubing kit.



The table below states the information of the Red tubing kit i.d. 0.13 mm (post-injection) for flow rates < 0.25 mL/min.

Label	Tubing	Length	Description
1A1	PEEK I.D. 0.25 mm	340 mm	Pump head to pump restrictor (left side)
1A2	PEEK I.D. 0.25 mm	340 mm	Pump head to pump restrictor (left side)
1B1	PEEK I.D. 0.25 mm	340 mm	Pump head to pump restrictor (right side)
1B2	PEEK I.D. 0.25 mm	340 mm	Pump head to pump restrictor (right side)
2A	PEEK I.D. 0.25 mm	115 mm	Pump restrictor (right side) to Pressure monitor (right side)
2B	PEEK I.D. 0.25 mm	115 mm	Pump restrictor (right side) to Pressure monitor (left side)
3	PEEK I.D. 0.25 mm	160 mm	Pressure monitor to Online filter
4	PEEK I.D. 0.25 mm	250 mm	Online filter to Injection valve V9M-J (port Pump)
5	PEEK I.D. 0.13 mm	520 mm	Injection valve V9M-J (port Col) to Column
(6)	M-M connector	N/A	Column to UV monitor U9-M
7	PEEK I.D. 0.13 mm	170 mm	UV monitor U9-M to Conductivity monitor C9M
8+9	PEEK I.D. 0.13 mm	175 mm	Conductivity monitor C9M to Outlet valve V9M-Os (port In)
Frac	PEEK I.D. 0.13 mm	400 mm	Outlet valve V9M-Os (Frac) to Fraction collector nozzle



[cytiva.com/chromatography](https://www.cytiva.com/chromatography)

Cytiva and the Drop logo are trademarks of Global Life Sciences IP Holdco LLC or an affiliate.

All other third-party trademarks are the property of their respective owners.

© 2020–2021 Cytiva

All goods and services are sold subject to the terms and conditions of sale of the supplying company operating within the Cytiva business. A copy of those terms and conditions is available on request. Contact your local Cytiva representative for the most current information.

For local office contact information, visit [cytiva.com/contact](https://www.cytiva.com/contact)

29380981 AC V:2 03/2021