

Injection Fill Port

Instructions

This instruction describes how to install and use the Injection Fill Port with ÄKTA™ injection valves (e.g., V9-Inj, V9H-Inj, V9-J, INV-907, and INV-917) for manual loading of sample into a sample loop. The Injection Fill Port has been optimized for use with glass syringes up to 2 mL. For larger volumes it is recommended to use a luer lock connection with a luer lock syringe.

Step	Action
1	Insert the Injection Fill Port nozzle into the injection valve port and lightly screw it on.
2	Insert the syringe needle into the Injection Fill Port. It is recommended to use a blunt ended 50 mm needle with a diameter of 0.70 mm.



Tip:

Leave a 2 mm gap between the Injection Fill Port and the syringe. This will make sure that any potential leak is detected.

- | | |
|---|------------------------------------------------------------------------------------------------------------------------------------------------------------|
| 3 | Tighten the Injection Fill Port until it forms a seal around the syringe needle.
Gently pull back the needle to make sure that it is securely fastened. |
|---|------------------------------------------------------------------------------------------------------------------------------------------------------------|

Step	Action
-------------	---------------

4	Inject your sample into the sample loop.
---	------------------------------------------

Note:

Leave the syringe in place until the sample is loaded onto the column to prevent back siphoning.

Page intentionally left blank



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

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

ÅKTA is a trademark of Global Life Sciences Solutions USA LLC or an affiliate doing business as Cytiva.

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)

29440098 AB V:9 04/2021