

42iQ Series pump

The purpose of this technical bulletin is to provide information related to the introduction of a new internal pump for the 42iQ series of NOx analyzers.

This new internal pump will be used in the models 42iQ, 42iQD, 42iQLS and 42iQHL (with bypass option).

The new pump is the same form, fit and function and meets all Thermo Scientific specifications and applicable certifications to be used in the above listed 42iQ series of NOx analyzers.

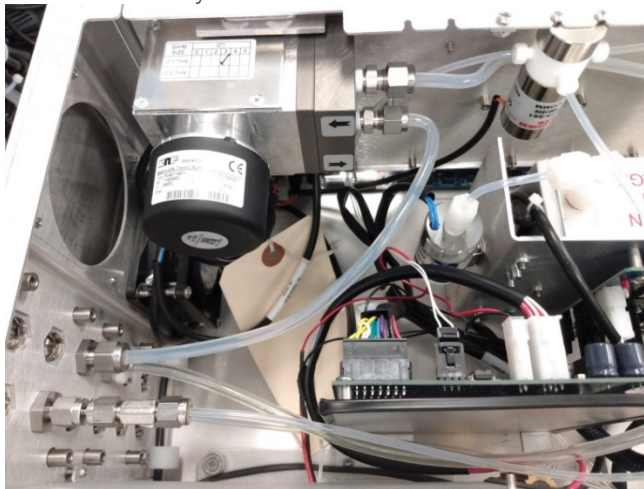


Image of new pump installed in 42iQ

There are two different ways of ordering the new pump,

- Customers replacing the existing double headed KNF DC pump use part number **120858-00 (Retro-fit kit)** as this will include the new pump, bulkhead fitting, Teflon tubing and installation instructions.
- If your 42iQ has shipped with or has been converted to this new pump and a replacement is needed, please order part number **120442-00 (Pump only)**.

New pump part numbers:

P/N: 120442-00 Pump only

P/N: 120858-00 Pump retrofit kit (Includes Pump, Bulkhead Fitting, PTFE Tubing and installation instruction)

P/N: 117901-00 Pump rebuild kit

USA

27 Forge Parkway

Franklin, MA 02038

Ph: (508) 520-0430

Toll Free: (866) 282-0430

orders.aqi@thermofisher.com

Find out more at [thermofisher.com](https://www.thermofisher.com)

© 2020 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.