

Recommended Holder Tightening for Prostak™ Tangential Flow Filtration (TFF) System Holders

The Prostak™ microfiltration modules utilize sealing gaskets between modules and endplates. It is designed such that it is not necessary to tighten to a specific torque, hand tightening is generally sufficient. For customers who prefer to use a torque wrench, it is recommended that they be uniformly tightened to 50–100 lbf-in.

If proper sealing does not occur, disassemble the holder carefully, verify that the gaskets are seated correctly and examine the Prostak™ modules for damage.

Avoid severely over torquing the filter modules as this may eventually crack the devices.



To place an order or receive technical assistance

In Europe, please call Customer Service:

France: 0825 045 645

Italy: 848 845 645

Switzerland: 0848 645 645

Germany: 069 86798021

Spain: 901 516 645 Option 1

United Kingdom: 0870 900 4645

For other countries across Europe, please call: +44 (0) 115 943 0840

Or visit: [MerckMillipore.com/offices](https://www.MerckMillipore.com/offices)

For Technical Service visit: [MerckMillipore.com/techservice](https://www.MerckMillipore.com/techservice)

[MerckMillipore.com](https://www.MerckMillipore.com)

The life science business of Merck operates as MilliporeSigma in the U.S. and Canada.

© 2018 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved. Merck, the vibrant M, Millipore and Prostak are trademarks of Merck KGaA, Darmstadt, Germany or its affiliates. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly accessible resources.

Lit. No. TB1004EN00 Ver. 2.0 2018-17332 11/2018