

Lab Close Kit

Lab Close Kit keeps your system in top condition when you're away

General Description

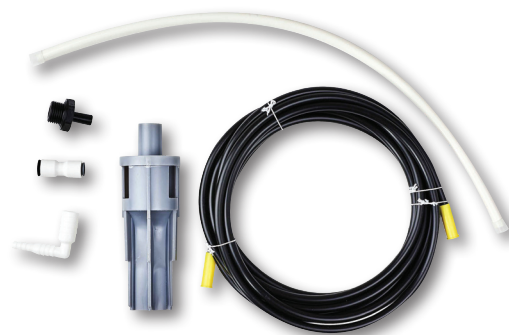
A kit that ensures your system will be automatically maintained during lengthy inactivity periods.

Key Benefits

- Minimizes water and energy usage during vacation periods or long inactivity periods
- Prevents contamination of water purification system and tank due to water stagnation
- Your system is up and running immediately when you come back

Ordering Information

Description	Cat. No.
Milli-Q® Integral; Elix® Advantage; Elix® Reference; and Milli-Q® Direct water purification systems	LABCLOSE1
RiOs™/Elix®/AFS® Essential water purification systems	included, setting to be activated



To place an order or receive technical assistance

In Europe, please call Customer Service:

France: 0825 045 645
Germany: 069 86798021
Italy: 848 845 645
Spain: 901 516 645 Option 1
Switzerland: 0848 645 645
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)
[merckgroup.com/life-science](https://www.merckgroup.com/life-science)

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, Milli-Q, AFS, Elix, RiOs, and the vibrant M 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.

MK_DS1251EN
2017 - 08107
1/2018