

Thermo Scientific Barnstead Pretreatment System Quick Reference Guide



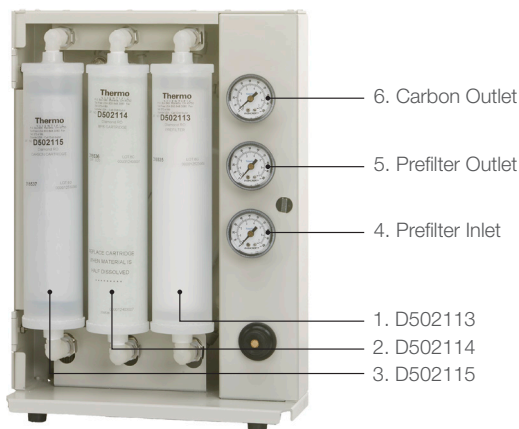
| Please refer to operation manual for complete instructions

Barnstead Pretreatment System removes feed water impurities that could harm reverse osmosis membranes. Use the table below for cartridge replacement frequency and pressure gauge monitoring.

Key#	Part #	Description	When to Replace
1	**D502113	1 micron prefilter - removes particles	<ul style="list-style-type: none"> • 6 months • Water flow to unit is reduced • Low pressure (see Key# 4 & 5)
2	**D502114	Hardness stabilizer - reduces water hardness	<ul style="list-style-type: none"> • Hardness stabilizer is half empty or 6 months, whichever comes first
3	**D502115	Carbon cartridge - removes chlorine	<ul style="list-style-type: none"> • 6 months • Chlorine is detected in concentrate water • Low pressure (see Key# 6) • Water flow at the outlet is reduced
Key#	It Monitors	It Measures	What to Monitor
4	Prefilter Inlet	Regulated water pressure before prefilter	<ul style="list-style-type: none"> • Pressure reading 0-15 psi, check feed water supply
5	Prefilter Outlet	Water pressure after the prefilter	<ul style="list-style-type: none"> • Pressure differential between Prefilter Inlet and Prefilter Outlet is >10 psi, change prefilter
6	Carbon Filter Outlet	Water pressure after the carbon filter	<ul style="list-style-type: none"> • Pressure differential between Prefilter Outlet and Carbon outlet is >15 psi or if pressure gauge is 0 psi.

**Included with new unit.

Note: The lifetime of consumables is a direct function of the quality of feed water and the daily volume of water that is used.



Find out more at thermofisher.com/purewater

©2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details. **50158356 0918**

ThermoFisher
SCIENTIFIC