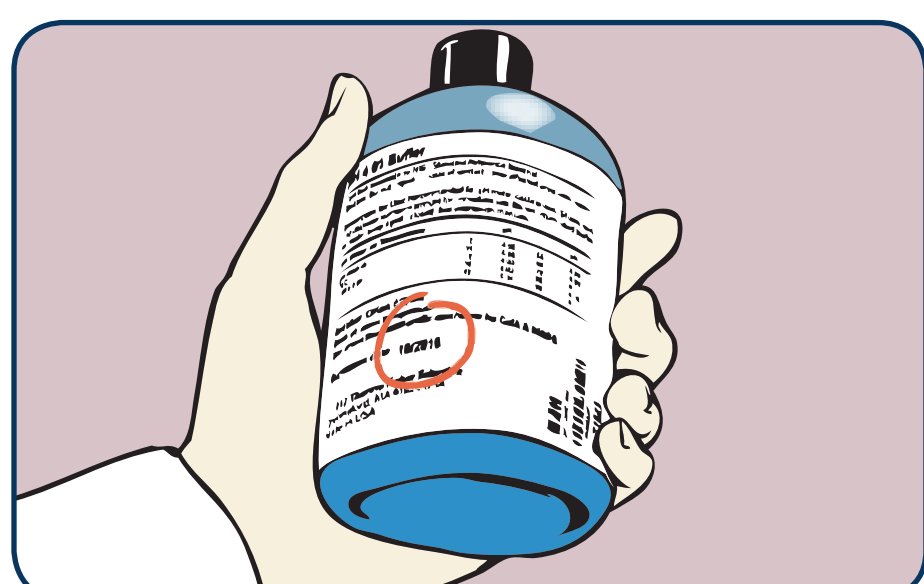


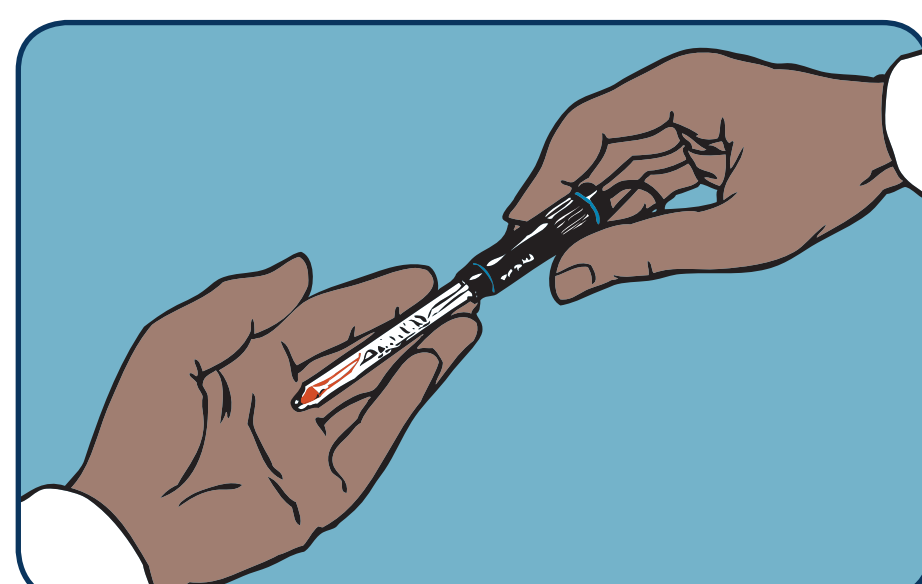


"Now then, when did you say was the last time you checked the expiration dates of these solutions?"

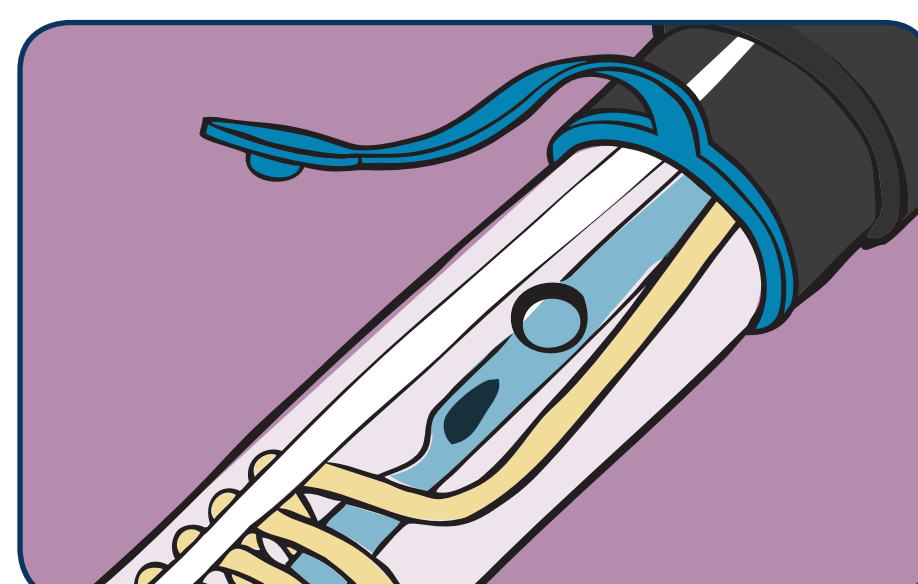
Good laboratory practices for the use, care and storage of Thermo Scientific Orion solutions and electrodes



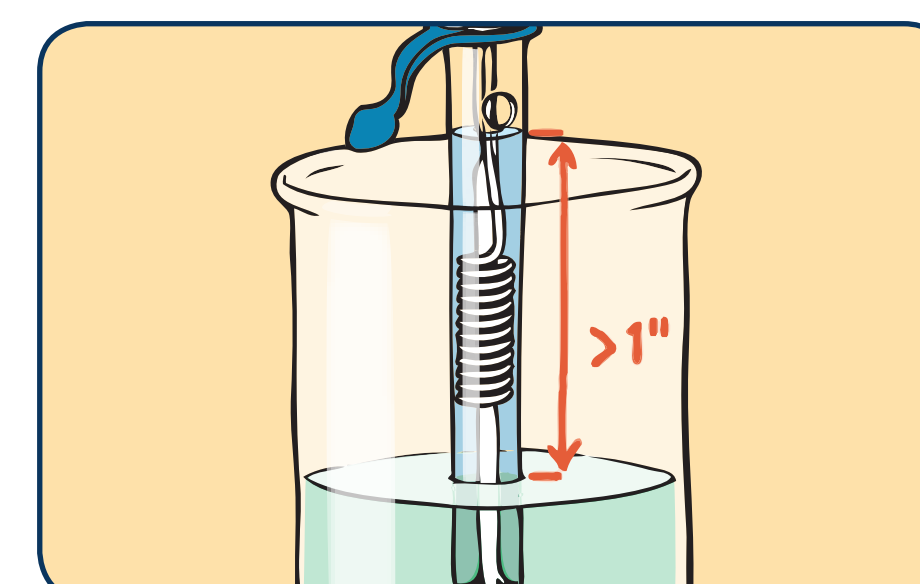
Always check the expiration date on the solution bottle prior to use. Unopened bottles of solution must be used within 2 years of the manufacture date. We recommend that opened bottles of solution be used within 1 - 3 months to avoid inaccurate measurements or changes in composition and calibration.



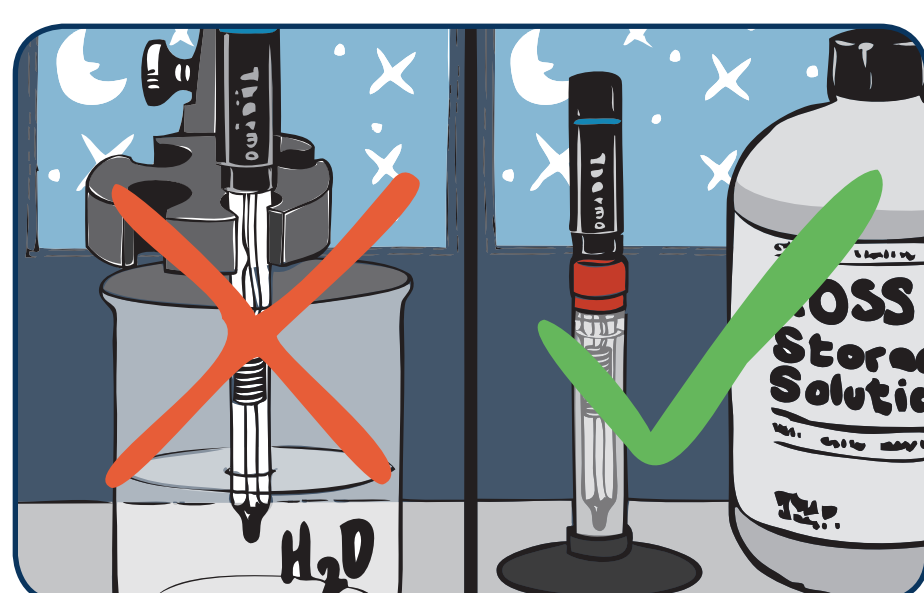
Periodically examine the surface of your electrode for crystallization near the fill hole, scratches or scuffs on the bulb surface, or discoloration of the filling solution. These may be signs that the electrode needs cleaning, replacement or new filling solution.



For best results with refillable electrodes, always keep the fill hole open during testing and closed for storage (use included fill hole cover for storage).



Fill the electrode to the bottom of the fill hole. Once filled, close the fill hole and gently shake the electrode to remove air bubbles. Open fill hole. When testing, the level of the fill solution should always be at least 1 inch (2.5 cm) above the level of sample being tested.



Deionized or distilled water is suitable for rinsing pH electrodes; however, neither should ever be used for long-term storage. When not in use, place the electrode in a sleeve half-filled with pH electrode storage solution. Never store an electrode dry.



To prevent carry-over contamination, rinse the electrode with deionized water. Never wipe the electrode bulb as a resulting static charge may interfere with the accuracy of the test. Either blot the bulb gently and briefly with a lint-free laboratory wipe, or gently shake excess water drops from the end.



Never put an unused portion of solution back into the bottle and do not reuse solution.



Store solutions between 60° and 85° F (15° and 30° C), out of direct sun and away from heat sources. Do not refrigerate.

Always use fresh solutions and properly maintained and functioning electrodes as good laboratory practice. Thermo Scientific™ Orion™ pH buffers and ISE standards have been specifically formulated to keep your Orion electrodes producing consistent, accurate results. Contact your authorized distributor to order new supplies.

thermoscientific.com/water

© 2014 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

PS1578-1001

Thermo
SCIENTIFIC

A Thermo Fisher Scientific Brand