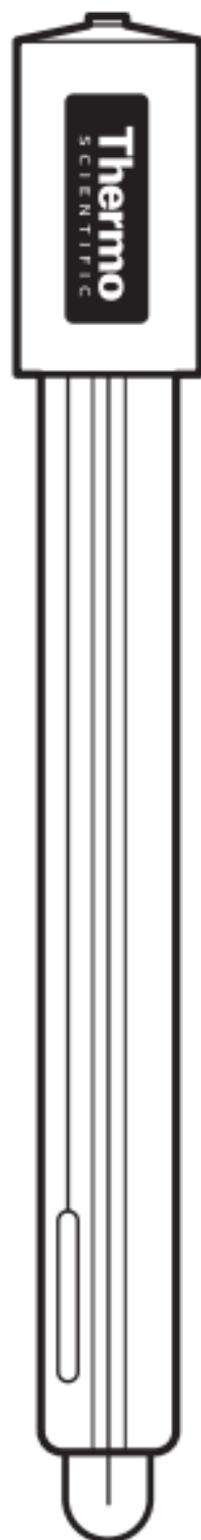


Glass pH Electrode 9142BN



- General purpose
- Glass refillable pH
- Combination electrode

9142BN

Lot Code:

OPERATION INSTRUCTIONS

Thank you for selecting this pH electrode.
pH range: 0-14. Dims (mm): 120 L x 12 dia.

BEFORE YOU BEGIN: Remove protective cap. Condition pH electrode, immerse it in fresh electrode storage solution or pH 4 or 7 buffer for 60 minutes before use.

CALIBRATION: Calibrate according to meter instruction. ALWAYS use fresh calibration solution. See recommended solution below.

DURING MEASUREMENT: To avoid contamination, rinse electrode with deionized water before and between measurement. Stir sample to ensure homogeneous mixture.

STORAGE: To store electrode, clean and dry electrode and replace protective cap.

ACCESSORIES: 900011 Filling Solution (5 x 60 mL); 910001 pH Storage Solution; 900020 pH Cleaning Solution Sample Kit; 910104 pH 4.01 buffer (475 mL); 910107 pH 7.00 buffer (475 mL); 910110 pH 10.01 buffer (475 mL).

WARRANTY: This electrode is warranted for a period of 6 months from the date of purchase. If replacement is necessary within the 6 month period and the electrode issue is not the result of abuse or misuse, please contact your local distributor.

RETURN OF ITEM: Authorization must be obtained before returning items for any reason. Note: Prices and design are subject to change without notice.

Made in USA

Distributed by Thermo Fisher Scientific

Environmental Instruments Water Analysis Instruments

166 Cummings Center, Beverly, MA 01915 USA
Toll Free: 1-800-225-1480 • Dom. Fax: 1-978-232-6015
Tel: 1-978-232-6000 • Int'l Fax: 978-232-6031

www.thermo.com/water