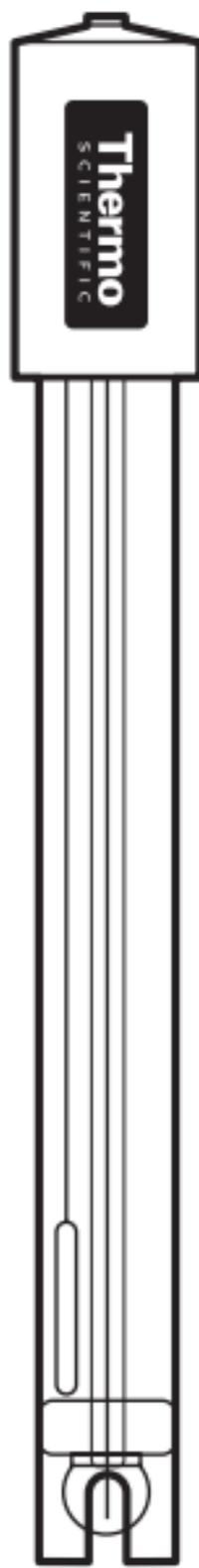


Gel pH Electrode 9146BN



- General purpose
- Non-refillable gel pH
- Combination electrode
- Rugged body design

9146BN

Lot Code:

Thermo
SCIENTIFIC

OPERATION INSTRUCTIONS

Thank you for selecting this pH electrode.
pH range: 0-14. Temp: 0-80 °C.
Dims (mm): 120 L x 12 dia.

BEFORE YOU BEGIN: Remove protective cap and place on bulb guard. To 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 solution to ensure homogeneous mixture.

STORAGE: To store electrode, clean and dry electrode. Remove bulb guard and replace with protective cap.

ACCESSORIES: 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 Singapore
Distributed by Thermo Fisher Scientific

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.thermoscientific.com/water