

Phosphatase Dilution Buffer I

Catalog # P21-09

Lot # S297-9

Product Description

Stock solution of buffer used for diluting phosphatase protein.

Formulation

Frozen solution of buffer containing: 25 mM HEPES pH 7.2, 50 mM NaCl, 2.5 mM EDTA, 65 ng/μl BSA.

Add to final concentration of 5 mM DTT prior to use. DTT Solution supplied in separate vial.

Sterility

Buffer passed through a 0.2μ filter membrane and packaged aseptically.

Storage and Stability

Store product at -20°C. For optimal storage, aliquot buffer into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

Phosphatase Dilution Buffer I

Catalog #

P21-09

Lot #

S297-9

Stability

1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot buffer into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped frozen.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.