

Catalog #	Aliquot Size
P23-09-01	1ml
P23-09-05	5ml
P23-09-20	20ml

Phosphatase Dilution Buffer III

Catalog # P23-09

Lot # U1873-4

Product Description

Solution of (1X) buffer used for diluting Phosphatase enzymes.

Formulation

Frozen solution of buffer containing: 20mM Tris-HCl, pH 7.5, 10mM MgCl₂, 0.02% Tween-20 and 65ng/μl BSA.

Add to final concentration of 0.4mM 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 III

Catalog #	P23-09
Lot #	U1873-4
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.