



Catalogue #	Aliquot Size
K22-09-01	1ml
K22-09-05	5ml
K22-09-20	20ml

Kinase Dilution Buffer II

Catalog # K22-09

Lot # S297-2

Product Description

Solution of buffer used for diluting kinase protein.

Formulation

Frozen solution of buffer containing: 5 mM MOPS, pH7.2, 2.5mM β -glycerol-phosphate, 4mM MgCl₂, 2.5mM MnCl₂, 1mM EGTA, 0.4mM EDTA

Add to final concentration of 0.05mM DTT (Catalog# D86-09B) prior to use. DTT Solution supplied in separate vial.

Sterility

Buffer passed through a 0.2 μ 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.

Kinase Dilution Buffer II

Catalog Number K22-09

Specific Lot Number S297-2

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.