



Catalog #	Aliquot Size
K25-09-01	1ml
K25-09-05	5ml
K25-09-20	20ml

Kinase Dilution Buffer V

Catalog # K25-09

Lot # L2039-8

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, 5mM MgCl₂, 0.4mM EDTA, 5% glycerol.

Add to final concentration of 50 μ M DTT (Catalog# D86-09) prior to use. DTT Solution supplied in separate vial.

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 V

Catalog #	K25-09
Lot #	L2039-8
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.