



| Catalog # | Aliquot Size |
|-----------|--------------|
| K26-09-01 | 1ml |
| K26-09-05 | 5ml |
| K26-09-20 | 20ml |

Kinase Dilution Buffer VI

Catalog # K26-09

Lot # L2039-9

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 $MgCl_2$, 0.4mM EDTA, 2.5mM $MnCl_2$, 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 $-20^{\circ}C$. 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 VI

| | |
|--------------------|---|
| Catalog # | K26-09 |
| Lot # | L2039-9 |
| Stability | 1yr at $-20^{\circ}C$ from date of shipment |
| Storage & Shipping | Store product at $-20^{\circ}C$. 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.