



Catalog #	Aliquot Size
K28-09-01	1ml
K28-09-05	5ml
K28-09-20	20ml

## **Kinase Dilution Buffer VIII**

**Catalog # K28-09**

Lot # L2039-11

### **Product Description**

Solution of buffer used for diluting kinase protein.

### **Formulation**

Frozen solution of buffer containing: 5 mM MOPS, pH7.2, 2.5mM  $\beta$ -glycerol-phosphate, 5mM  $MgCl_2$ , 0.4mM EDTA, 2.5mM  $MnCl_2$ , 50ng/ $\mu$ l BSA, 5% glycerol.

Add to final concentration of 50 $\mu$ 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 VIII**

Catalog #	K28-09
Lot #	L2039-11
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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

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