



Catalog #	Aliquot Size
K23-09-01	1ml
K23-09-05	5ml
K23-09-20	20ml

## Kinase Dilution Buffer III

Catalog # K23-09

Lot # S297-3

### 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  $MgCl_2$ , 1mM EGTA, 0.4mM EDTA, 50ng/ $\mu$ l BSA.

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 $\mu$  filter membrane and packaged aseptically.

### 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 III

Catalog Number K23-09  
Specific Lot Number S297-3

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.**