

## Kinase Dilution Buffer V

**Catalog # K25-09**

Lot # S297-5

### 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<sub>2</sub>, 1mM EGTA, 0.4mM EDTA, 0.05mM DTT, 5% glycerol.

### Sterility

Buffer passed through a 0.2 $\mu$  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 V

Catalog Number K25-09

Specific Lot Number S297-5

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](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.**