

## Kinase Assay Buffer II

**Catalog # K02-09**

Lot # S285-2

### Product Description

Stock solution of buffer used for assaying kinase protein.

### Formulation

Frozen solution of buffer containing: 25 mM MOPS, pH7.2, 12.5mM  $\beta$ -glycerol-phosphate, 20mM MgCl<sub>2</sub>, 12.5mM MnCl<sub>2</sub>, 5mM EGTA, 2mM EDTA.

Add to final concentration of 0.25mM 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°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 Assay Buffer II

Catalog Number K02-09

Specific Lot Number S285-2

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.**