



Catalog #

Aliquot Size

K20-09-01  
K20-09-05  
K20-09-20

1ml  
5ml  
20ml

## Kinase Dilution Buffer X

**Catalog # K20-09**

Lot # M2839-6

### Product Description

The working solution used for diluting kinase proteins.

### Formulation

Frozen solution of buffer containing: 40mM Tris-HCl, pH7.4, 20mM MgCl<sub>2</sub>, 2.5mM MnCl<sub>2</sub>, and 0.1mg/ml BSA.

Add to final concentration of 50µM DTT (Catalog# D86-09) to the aliquot prior to use. DTT Solution supplied in separate vial.

### Sterility

Buffer passed through a 0.2µ 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 X

Catalog #

K20-09

Lot #

M2839-6

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.**