

## Phosphatase Assay Buffer I

**Catalog # P01-09**

Lot # S285-3

### Product Description

Stock solution of buffer used for assaying phosphatase protein.

### Formulation

Frozen solution of buffer containing: 125 mM HEPES pH 7.2, 250 mM NaCl, 12.5 mM EDTA.

### Sterility

Buffer passed through a 0.2 $\mu$  filter membrane and packaged aseptically.

### Storage and Stability

Store product at  $-20^{\circ}\text{C}$ . 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.

## Phosphatase Assay Buffer I

Catalog Number P01-09

Specific Lot Number S285-3

Stability 1yr At  $-20^{\circ}\text{C}$  from date of shipment  
Storage & Shipping Store product at  $-20^{\circ}\text{C}$ . 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.**