

## Phosphatase Dilution Buffer II

**Catalog # P22-09**

Lot # S297-10

### Product Description

Stock solution of buffer used for diluting phosphatase protein.

### Formulation

Frozen solution of buffer containing: 50 mM Imidazole, pH 7.2, 0.2% 2-mercaptoethanol, 65ng/ $\mu$ l BSA.

### 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 Dilution Buffer II

Catalog Number P22-09

Specific Lot Number S297-10

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.**