

Catalog #	Aliquot Size
P03-09-01	1ml
P03-09-05	5ml
P03-09-20	20ml

## Phosphatase Assay Buffer III

**Catalog # P03-09**

Lot # U1873-3

### Product Description

5X stock solution of the buffer used for assaying phosphatase proteins.

### Formulation

Frozen solution of buffer containing: 100mM Tris-HCl pH 7.5, 50mM MgCl<sub>2</sub> and 0.1% Tween-20.

Add to final concentration of 2mM DTT (Catalog# D86-09B) and 300ng/μl BSA 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.

## Phosphatase Assay Buffer III

Catalog	P03-09
Specific Lot	U1873-3
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.**