

## Methyltransferase Dilution Buffer II

**Catalog # M22-09**

Lot # O1082-8

### Product Description

Solution of buffer used for diluting enzyme protein methyltransferases.

### Formulation

Frozen solution of buffer containing: 50mM Tris-HCl, pH 8.0 and 10ng/μl BSA.

Add to final concentration of 0.5mM DTT (Catalog# D86-09B) 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.

## Methyltransferase Dilution Buffer II

Catalog Number M22-09

Specific Lot Number O1082-8

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.**