

Catalogue #	Aliquot Size
M03-09-01	1ml
M03-09-05	5ml
M03-09-20	20ml

## **Methyltransferase Assay Buffer III**

**Catalog # M03-09**

Lot # O1035-2

### **Product Description**

5X stock solution of the buffer used for assaying DNA methyltransferases.

### **Formulation**

Frozen solution of buffer containing: 250mM Tris-HCl, pH 7.5, 5mM EDTA, 25% Glycerol, 50ng/μl BSA.

Add to final concentration of 2.5mM 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 Assay Buffer III**

Catalog Number	M03-09
Specific Lot Number	O1035-2
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.**