

Catalogue #	Aliquot Size
M23-09-01	1ml
M23-09-05	5ml
M23-09-20	20ml

Methyltransferase Dilution Buffer III

Catalog # M23-09

Lot # O1035-3

Product Description

Solution of (1X) buffer used for diluting enzyme DNA methyltransferases.

Formulation

Frozen solution of buffer containing: 50mM Tris-HCl, pH 7.5, 1mM EDTA, 5% Glycerol and 10ng/ μ l 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 0.2 μ 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 III

Catalog Number M23-09

Specific Lot Number O1035-3

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
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.