

Catalogue #	Aliquot Size
M21-09-01	1ml
M21-09-05	5ml
M21-09-20	20ml

Methyltransferase Dilution Buffer I

Catalog # M21-09

Lot # O1082-7

Product Description

Solution of buffer used for diluting enzyme protein methyltransferases.

Formulation

Frozen solution of buffer containing: 50mM Tris-HCl, pH 9.0 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 I

Catalog Number M21-09

Specific Lot Number O1082-7

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.