

Catalogue #	Aliquot Size
M02-09-01	1ml
M02-09-05	5ml
M02-09-20	20ml

Methyltransferase Assay Buffer II

Catalog # M02-09

Lot # O1082-6

Product Description

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

Formulation

Frozen solution of buffer containing: 250mM Tris-HCl, pH 8.0 and 50ng/ μ l 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 0.2 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 Assay Buffer II

Catalog Number M02-09

Specific Lot Number O1082-6

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.