



Catalog #	Aliquot Size
K03-09-01	1ml
K03-09-05	5ml
K03-09-20	20ml

Kinase Assay Buffer III

Catalog # K03-09

Lot # L2183-9

Product Description

The 5x stock solution used for assaying kinase proteins.

Formulation

Frozen solution of buffer containing: 200mM Tris-HCl, pH7.4, 100mM MgCl₂ and 0.5mg/ml BSA.

Add to final concentration of 0.25mM DTT (Catalog# D86-09) to the aliquot prior to use. 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.

Kinase Assay Buffer III

Catalog #	K03-09
Lot #	L2183-9
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.