

Lipid Kinase Buffer

Catalog # L01-09

Lot # L2111-15

Product Description

The 5x stock solution of buffer used for assaying lipid kinase proteins.

Formulation

Frozen stock solution of buffer containing: 250mM HEPES, pH7.5, 250mM NaCl, 10mM MgCl₂, 2.5mM EGTA, 125µg/mL BSA and 0.2% Triton X-100.

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 0.2µ 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.

Lipid Kinase Buffer

Catalog #	L01-09
Lot #	L2111-15
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.