

AMP Solution

Catalog # A46-09

Lot # X406-4

Product Description

0.5mM AMP solution used for AMPK family kinase assays.

Formulation

Frozen solution of buffer containing: 25mM MOPS pH 7.2, and 0.5mM AMP.

Sterility

Solution 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.

AMP Solution

Catalog Number A46-09

Specific Lot Number X406-4

Concentration 0.5 mM

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.