



Catalog #

Aliquot Size

A50-09-200

200ul

ATP Stock Solution

Catalog # A50-09

Lot # A1328-8

Product Description

10 mM ATP Stock Solution used for preparing hot ATP assay cocktail.

Formulation

Frozen solution containing: 10 mM ATP, 25 mM MOPS, pH 7.2, 12.5mM β -glycerol-phosphate, 25mM $MgCl_2$, 2mM EDTA and 0.25mM DTT.

Sterility

Buffer passed through a 0.2 μ filter membrane and packaged aseptically.

Storage and Stability

Store product at $-20^{\circ}C$. 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.

ATP Stock Solution

Catalog # A50-09

Lot # A1328-8

Concentration 10mM

Stability 1yr at $-20^{\circ}C$ from date of shipment

Storage & Shipping Store product at $-20^{\circ}C$. 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.