

GTP Solution

Catalog # G50-09

Lot # W175-5

Product Description

500 μ l volume of 5 μ M GTP solution for different enzyme assays.

Formulation

Frozen solution of buffer containing: 25mM MOPS pH 7.2, and 5 μ M GTP.

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.

GTP Solution

Catalog #	G50-09
Volume	500 μ l
Lot #	W175-5
Concentration	5 μ M
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.