

cGMP Solution

Catalog # G47-09

Lot # W175-4

Product Description

500µl volume of 100µM cGMP solution for different enzyme assays.

Formulation

Frozen solution of buffer containing: 25mM MOPS pH 7.2, and 100µM cGMP.

Sterility

Solution 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.

cGMP Solution

Catalog Number G47-09

Volume 500 µl

Specific Lot Number W175-4

Concentration 100µM

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.