

DLG (Dilauroyl-sn-glycerol)

Catalog # D430-59

Lot # M2900-7

Product Description

The substrate solution used for assaying Diacylglycerol kinases.

Molecular Weight

The molecular weight of 1,2-Dilauroyl-sn-glycerol (DLG) is: 456.70.

Formulation

Frozen substrate solution containing: 10 mM of 1,2-Dilauroyl-sn-glycerol (DLG) in DMSO.

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.

DLG (Dilauroyl-sn-glycerol)

Catalog #	D430-59
Lot #	M2900-7
Concentration	10 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.