



Catalog #

Aliquot Size

P422-59-500
P422-59-1000

0.5 ml
1.0 ml

PI(4)P

Catalog # P422-59

Lot # L2111-7

Product Description

The substrate solution used for assaying lipid kinases.

Molecular Weight

The molecular weight of Phosphatidylinositol 4-phosphate [PI(4)P or PtdIns4P] is: 620.32.

Formulation

Frozen substrate solution containing: 1mM of Phosphatidylinositol 4-phosphate [PI(4)P or PtdIns4P] in 1x Lipid Dilution Buffer (Catalog# L21-09).

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.

PI(4)P

Catalog #

P422-59

Lot #

L2111-7

Concentration

1 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.