

PI(4,5)P2:PS

Catalog # P429-59

Lot # L2111-14

Product Description

The substrate solution used for assaying lipid kinases.

Molecular Weight

The molecular weight of Phosphatidylserine (PS) is: 812.05;
The molecular weight of Phosphatidylinositol 4,5-bi-phosphate [PI(4,5)P2 or PtdIns(4,5)P] is: 856.48.

Formulation

Frozen substrate solution containing: 0.125mM of Phosphatidylinositol 4,5-biphosphate [PI(4,5)P2 or PtdIns(4,5)P] and 1mM of Phosphatidylserine (PS) 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(4,5)P2:PS

Catalog #

P429-59

Lot #

L2111-14

Stability

Storage & Shipping

1yr at -20°C from date of shipment
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.