

PRODUCT INFORMATION



PtdIns-(4,5)-P₂-fluorescein (triethylammonium salt)

Item No. 10010388

Formal Name: 1-(1-octadecanoyl-fluorescein-2R-octadecanoylphosphatidyl)inositol-4,5-bisphosphate, pentatriethylammonium salt

Synonyms: DOPI-4,5-P₂-fluorescein, Phosphatidylinositol-4,5-diphosphate C-8-fluorescein, PI(4,5)P₂-fluorescein, PIP2[4',5']-fluorescein

MF: C₄₆H₅₅NO₂₅P₃ • 5C₆H₁₆N

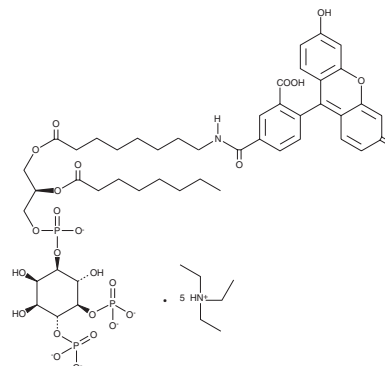
FW: 1,625.9

Purity: ≥95%

Supplied as: A lyophilized powder

Storage: -20°C

Stability: ≥1 year



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

PtdIns-(4,5)-P₂-fluorescein (triethylammonium salt) is supplied as a lyophilized powder. The solubility of PtdIns-(4,5)-P₂-fluorescein (triethylammonium salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

The phosphatidylinositol (PtdIns) phosphates represent a small percentage of total membrane phospholipids. However, they play a critical role in the generation and transmission of cellular signals.^{1,2} PtdIns-(4,5)-P₂-fluorescein is a fluorescent probe for any protein with a high-affinity binding interaction with inositol-(4,5)-triphosphate phospholipids, such as PI3K, PTEN, or PH-domain-containing proteins.

References

1. Exton, J.H. Regulation of phosphoinositide phospholipases by hormones, neurotransmitters, and other agonists linked to G proteins. *Annu. Rev. Pharmacol. Toxicol.* **36**, 481-509 (1996).
2. Majerus, P.W. Inositol phosphate biochemistry. *Annu. Rev. Biochem.* **61**, 225-250 (1992).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 09/21/2023

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM