

PRODUCT INFORMATION



Sulforhodamine B-Phosphatidylethanolamine (ammonium salt)

Item No. 45093

CAS Registry No.: 383906-59-0

Synonym: Sulforhodamine B-PE

MF: $C_{66}H_{103}N_3O_{14}PS_2 \cdot NH_4$

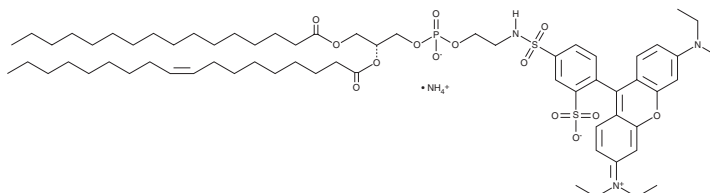
FW: 1,275.7

Purity: $\geq 95\%$ (mixture)

Supplied as: A 1 mg/ml solution in chloroform

Storage: $-20^\circ C$

Stability: ≥ 4 years



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

Description

Sulforhodamine B-phosphatidylethanolamine is a phosphatidylethanolamine isolated from chicken egg that has been labeled with an N-linked fluorescent sulforhodamine B (SRB; Item No. 31539) group. Sulforhodamine B-phosphatidylethanolamine has been used in membrane fusion assays and in the study of *in vitro* and *in vivo* liposome delivery and localization.¹⁻³ Sulforhodamine B displays excitation/emission maxima of 560/583 nm, respectively. This product contains phosphatidylethanolamine molecular species isolated from chicken egg with primarily C16:0, C18:0, C18:1, C18:2, and C20:4 fatty acyl chains.

References

1. Loura, L.M.S. and Prieto, M. FRET in membrane biophysics: An overview. *Front. Physiol.* **2**, 82 (2011).
2. Lymberopoulos, A., Demopoulou, C., Kyriazi, M., *et al.* Liposome percutaneous penetration in vivo. *Toxicol. Res. Appl.* **1**, 1-6 (2017).
3. Hazemoto, N., Harada, M., Komatsubara, N., *et al.* pH-sensitive liposomes composed of phosphatidylethanolamine and fatty acid. *Chem. Pharm. Bull. (Tokyo)* **38(3)**, 748-751 (1990).

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, 07/06/2026

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