## **Rhodamine-DHPE**

Rhodamine-DHPE is a rhodamine-labeled glycerophosphoethanolamine lipid that has been used in membrane fusion assays.



## **Product attributes**

CAS number	126111-99-7
Probe cellular localization	Membrane/cell surface, Membrane/vesicular
Colors	Red
Excitation/Emission	560/581 nm
Conjugation	Rhodamine

Call us: 800-304-5357 Email: btinfo@biotium.com

## **Product Description**

Rhodamine-DHPE (N-(Lissamine rhodamine B

sulfonyl)-1,2-dihexadecanoyl-sn-glycero-3-phosphoethanolamine, triethylammonium salt) is a rhodamine-labeled glycerophosphoethanolamine lipid. DHPE conjugates of rhodamine dyes have been used as fluorescence energy acceptors in combination with NBDPE as the donor in membrane fusion assays using the principle of fluorescence energy transfer (FRET) (1-3). These probes have also been used for following membrane trafficking during endocytosis (4,5).

- $\lambda_{Ex}/\lambda_{Em}$  (MeOH) = 560/581 nm
- Dark red solid soluble in chloroform
- Store at -20°C and protect from light
- $\bullet \ C_{70}H_{117}N_4O_{14}PS_2$
- MW: 1333.80
- [126111-99-7]

$$(\text{CH}_3\text{CH}_2\text{C})_2\text{N} \\ (\text{CH}_3\text{CH}_2)_3\text{N}\text{H} \\ O \\ \text{H}_3\text{C}(\text{H}_2\text{C})_{14} \\ -\text{C} \\ -\text{OCH}_2 \\ \text{H}_3\text{C}(\text{H}_2\text{C})_{14} \\ -\text{C} \\ -\text{OCH}_2 \\ \text{O} \\ \text{CH}_2\text{O} \\ -\text{P} \\ -\text{OCH}_2\text{CH}_2\text{NH} \\ \text{O} \\ \end{array}$$

## References

- 1. Biochemistry 24, 6390 (1985).
- 2. Biochemistry 21, 1720 (1982).
- 3. Biochemistry 20,

4093 (1981).

- 4. Biochim Biophys Acta 1103, 185 (1992).
- 5. Eur J Cell Biol 53, 173 (1990).

This datasheet was generated on November 3, 2025 at 06:03:32 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/rhodamine-triethylammonium-salt/">https://biotium.com/product/rhodamine-triethylammonium-salt/</a>