

Sphingosine-tetramethylrhodamine

Catalog number: S-100T

Molecular Formula: C₄₃H₅₉N₄O₆

MW: 727.95

CAS#: n/a

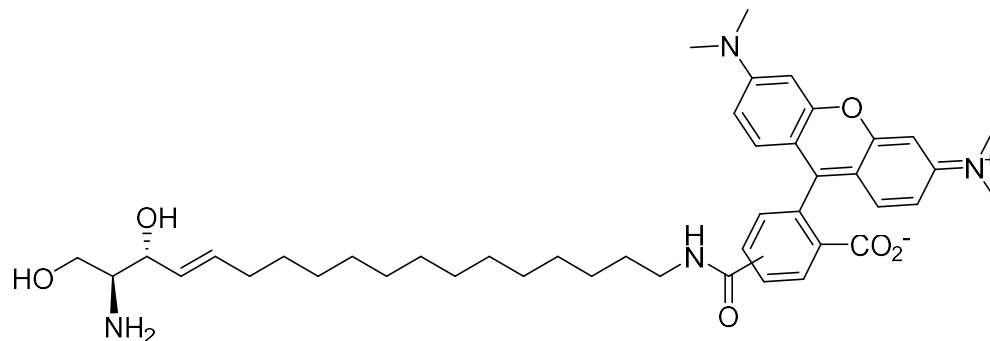
Excitation/Emission: 555/580 nm

Solubility: MeOH, >1 mg/ml. Stock solutions can be further diluted into aqueous buffer.

Storage and Handling: Sphingosine-tetramethylrhodamine is relatively stable at room temperature as a solid. It is best stored as a lyophilized solid at -20°C. Solutions/emulsions of sphingosines should be kept at -20°C. Storage in basic buffers (pH > 9.0) or acidic buffers (pH < 4.0) may cause decomposition. Protect from light.

Background: Sphingosine is formed through ceramidase catalyzed hydrolysis of ceramide. It is a potent inhibitor of protein kinase C (IC₅₀ = 1-3 μM). Sphingosine is phosphorylated by sphingosine kinase to the bioactive lipid sphingosine 1-phosphate.

References: 1) C.R. Gault, *et al.* (2010) "An overview of sphingolipid metabolism: from synthesis to breakdown" *Adv. Exp. Med. Biol.* 688: 1-23.
2) Y. A. Hannum & L.M. Obeid (2017) "Sphingolipids and their metabolism in physiology and disease" *Nat. Rev. Mol. Cell Biol.* 19:175-191.
3) M. Rekhter & S. Karathanasis (2006) "Sphingolipids in atherosclerosis: a metabolic underpinning of vascular disease" *Future Lipidology* 1(5): 204-614.
4) Lloyd-Evans, E., A. J. Morgan, *et al.* (2008). "Niemann-Pick disease type C1 is a sphingosine storage disease that causes deregulation of lysosomal calcium." *Nat Med* 14(11): 1247.



Hazardous Properties and Cautions: The toxicological and pharmacological properties of this compound are not fully known. For further information see the MSDS on request. This product is manufactured and shipped only in small quantities, intended for research and development in a laboratory utilizing prudent procedures for handling chemicals of unknown toxicity, under the supervision of persons technically qualified to evaluate potential risks and authorized to enforce appropriate health and safety measures. As with all research chemicals, precautions should be taken to avoid unnecessary exposures or risks.

Warranty and Disclaimer: Echelon warrants the product conforms to the specifications stated herein. In the event of nonconformity, Echelon will replace products or refund purchase price, at its sole option, and Echelon shall not be responsible for any other loss or damage, whether known or foreseeable to Echelon. No other warranties apply, express or implied, including but not limited to warranty of fitness for any purpose or implied warranty of merchantability. Purchaser is solely responsible for all consequences of its use of the product and Echelon assumes no responsibility therefore, including success of purchaser's research and development, or health or safety of any uses of the product.

Technical Data Sheet, Rev 6, 5/13/24– **For research use only.** Not intended for diagnostic or therapeutic use.