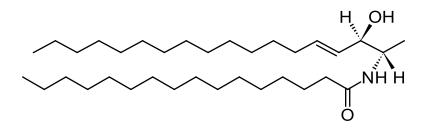
TECHNICAL DATA SHEET

N-palmitoyl-D-erythro-sphingosine (d18:1/16:0)

Catalog Number	860456	Physical state	Powder
Purity	> 99%	Transition temp.	No data
CAS		СМС	No data
Synonyms	N-C16-deoxy-sphingosine; N-hexa- decanoyl-D-erythro-sphing-4-enine (d18:1/16:0)	pK _a	No data
Molec. Formula	C ₃₄ H ₆₇ NO ₂	TLC mobile phase	C:M*, 95:5, v/v
MW	521.901	Exact Mass	521.517
Percent composition	C 78.25% H 12.94% N 2.68% O 6.13%		
Stability	Store in <-20°C freezer for up to one year		
Solubility	Soluble in DMSO, Methanol, Ethanol:Water (95:5) at 40°C + Sonication. Note: solidifies at room temperature in all solvents (i.e. keep warm with sonication). Concentration should be 5 mg/mL in DMSO; 10 mg/mL in methanol or ethanol:water (95:5).		
Web link	860456		

*C, chloroform; M, methanol



Description:

Commonly referred to as 1-deoxydihydroceramide (1-deoxyDHCer), this product is the N-acylated form of 1-deoxysphinganine, a potent inhibitor of sphingolipid metabolism. The biological activity of 1-deoxyDHCer is not clearly understood at this time.

References:

Zitomer, NC *et al* (2009) Ceramide synthase inhibition by fumosin B1 causes accumulation of 1-deoxysphinganine. J Biol Chem. 284(8):4786-95
Pruett, ST *et al* (2008). Thematic Review Series: Sphingolipids. Biodiversity of sphingoid bases ("sphingosines") and related amino alcohols. J Lipid Res 49:1621-39

Related products: Sphingolipids

MSDS: Available at www.avantilipids.com for Product Number 860456

Avanti Polar Lipids, Inc., 700 Industrial Park Drive, Alabaster, AL 35007-9105, U.S.A. Phone •(800) 227-0651 •(205) 663-2494 •Fax (800) 229-1004 •(205) 663-0756 Email •orders@avantilipids.com •Technical Questions: technical@avantilipids.com •Inquiries: info@avantilipids.com •Analytical: analytical@avantilipids.com

