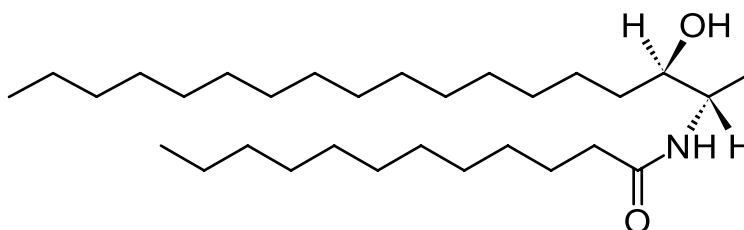


TECHNICAL DATA SHEET

N-lauroyl-1-deoxy-D-erythro-sphinganine (d18:0/12:0)

Catalog Number	860460	Physical state	Powder
Purity	> 99%	Transition temp.	No data
CAS		CMC	No data
Synonyms	N-C12-deoxy-sphinganine; N-dodecanoyl-1-deoxy-D-erythro-sphinganine (d18:0/12:0)	pK_a	No data
Molec. Formula	C ₃₀ H ₆₁ NO ₂	TLC mobile phase	C:M*, 95:5, v/v
MW	467.811	Exact Mass	467.470
Percent composition	C 77.02% H 13.14% N 2.99% O 6.84%		
Stability	Store in <-20°C freezer for up to one year		
Solubility	Soluble in DMSO at 40°C + Sonication. Note: solidifies at room temperature in DMSO (i.e. keep warm with sonication). Soluble in methanol and ethanol:water, 95:5, at 40°C. Concentration should be 10 mg/mL.		
Web link	860460		

*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 860460