

PRODUCT DATA SHEET

3-keto-C8-Dihydroshingosine·HCl

Catalog number: 1892

Common names: 1-Hydroxy-2-amino-3-keto-octane·HCl

Source: synthetic

Solubility: chloroform, methanol, ethanol, DI water

CAS number: 1824382-78-6

Molecular Formula: C₈H₁₇NO₂·HCl

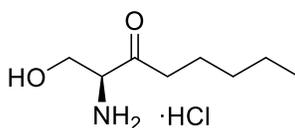
Molecular Weight: 159

Storage: -20°C

Purity: TLC: 98%

TLC System: chloroform/methanol (80:20)

Appearance: solid



Application Notes:

3-keto-Dihydroshingosine is a vital intermediate in the biosynthesis of ceramides.^{1,2} This lyso-sphingolipid is formed by the condensation of L-serine and palmitoyl-CoA by the serine palmitoyl transferase enzyme.³ It is then reduced to dihydroshingosine, converted to ceramide, and eventually synthesized into many types of sphingolipids. C18-keto-dihydroshingosine is the major sphingolipid precursor in the early and intermediate stages of cell life with C20-keto-dihydroshingosine as a minor component. Towards the end of the cell's life the ratio of C18 to C20 keto-dihydroshingosine becomes more equal. It seems to be a critical regulating step in the availability of sphingolipids in cells. Vitamin K deficiency results in the inactivation of the serine palmitoyl transferase enzyme causing a resultant shortage of sphingolipids. Short-chain analogs of 3-keto-dihydroshingosine have different physical properties from the long-chain 3-keto-dihydroshingosines. Short-chain bases are considerably less hydrophobic which could significantly change the process of signal transduction.

Selected References:

1. N. Bartke and Y. Hannun "Bioactive sphingolipids: metabolism and function" *Journal of Lipid Research*, Vol. 50 pp. S91-S96, 2009
2. G. Jenkins and Y. Hannun "Role for *de Novo* Sphingoid Base Biosynthesis in the Heat-induced Transient Cell Cycle Arrest of *Saccharomyces cerevisiae*" *Journal of Biological Chemistry*, Vol. 276 pp. 8574-8581, 2001
3. A. Batheja et al. "Characterization of Serine Palmitoyltransferase in Normal Human Tissues" *Journal of Histochemistry and Cytochemistry*, Vol. 51 pp. 687-696, 2003

This product is to be used for research only. It is not intended for drug or diagnostic use, human consumption or to be used in food or food additives. Matreya assumes no liability for any use of this product by the end user. We believe the information, offered in good faith, is accurate.

DS1892 Rev. #5
November 2, 2021