

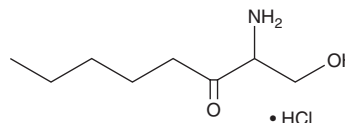
# PRODUCT INFORMATION



## 3-keto Sphinganine (d8:0) (hydrochloride)

Item No. 22824

**CAS Registry No.:** 1824382-78-6  
**Formal Name:** 2-amino-1-hydroxy-3-octanone, monohydrochloride  
**Synonym:** 3-keto-C8-Dihydrosphingosine  
**MF:** C<sub>8</sub>H<sub>17</sub>NO<sub>2</sub> • HCl  
**FW:** 195.7  
**Purity:** ≥98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

3-keto Sphinganine (d8:0) (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the 3-keto Sphinganine (d8:0) (hydrochloride) in the solvent of choice, which should be purged with an inert gas. 3-keto Sphinganine (d8:0) (hydrochloride) is soluble in organic solvents such as ethanol, methanol, and chloroform. 3-keto Sphinganine (d8:0) (hydrochloride) is also soluble in DI water. We do not recommend storing the aqueous solution for more than one day.

### Description

3-keto Sphinganine (d8:0) is a short-chain analog of the typical C18 chain-length 3-keto sphinganine (d18:0) (Item No. 24380). 3-keto Sphinganine (d18:0) is a lyso-sphingolipid formed by the condensation of L-serine and palmitoyl-CoA by serine palmitoyl transferase (SPT). Vitamin K deficiency inactivates SPT resulting in a decrease in the production of 3-keto sphinganine and other sphingolipids.<sup>1</sup>

### References

1. Batheja, A.D., Uhlinger, D.J., Carton, J.M., *et al.* Characterization of serine palmitoyltransferase in normal human tissues. *J. Histochem. Cytochem.* **51(5)**, 687-696 (2003).

**WARNING**  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

**SAFETY DATA**  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

**WARRANTY AND LIMITATION OF REMEDY**  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/16/2022

### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM