

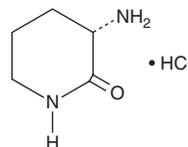
PRODUCT INFORMATION



L-Ornithine lactam (hydrochloride)

Item No. 31683

CAS Registry No.: 42538-31-8
Formal Name: 3S-amino-2-piperidinone, monohydrochloride
MF: $C_5H_{10}N_2O \cdot HCl$
FW: 150.6
Purity: $\geq 95\%$
Supplied as: A crystalline solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



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

Laboratory Procedures

L-Ornithine lactam (hydrochloride) is supplied as a crystalline solid. Aqueous solutions of L-ornithine lactam (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of L-ornithine lactam (hydrochloride) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

L-Ornithine lactam is a synthetic intermediate.^{1,2} It has been used in the synthesis of dipeptidyl peptidase 4 (DPP-4) inhibitors, as well as signaling peptides.

References

1. Abe, M., Akiyama, T., Umezawa, Y., *et al.* Synthesis and biological activity of sulphostin analogues, novel dipeptidyl peptidase IV inhibitors. *Bioorg. Med. Chem.* **13(3)**, 785-797 (2005).
2. Barnett, R., Raszkowski, D., Winckler, T., *et al.* Versatile synthesis of the signaling peptide glorin. *Beilstein J. Org. Chem.* **13**, 247-250 (2017).

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, 11/14/2022

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