

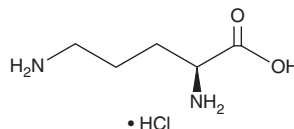
PRODUCT INFORMATION



L-Ornithine (hydrochloride)

Item No. 30779

CAS Registry No.: 3184-13-2
Formal Name: L-ornithine, monohydrochloride
MF: $C_5H_{12}N_2O_2 \cdot HCl$
FW: 168.6
Purity: $\geq 95\%$
Supplied as: A 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 (hydrochloride) is supplied as a solid. Aqueous solutions of L-ornithine (hydrochloride) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of L-ornithine (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 is an active metabolite of L-arginine (Item No. 23703) and a non-proteinogenic amino acid.¹ L-Ornithine is formed from arginine by arginase I during the final step of the urea cycle and acts as an alternate precursor to L-glutamate in the biosynthesis of proline. Exogenous administration of L-ornithine (111 mg/kg) delays the onset of ethanol-induced ataxia and decreases ethanol-induced sleeping time in rats.²

References

1. Häberle, J., Boddaert, N., Burlina, A., *et al.* Suggested guidelines for the diagnosis and management of urea cycle disorders. *Orphanet J. Rare Dis.* 7, 32 (2012).
2. Breglia, R.J., Ward, C.O., and Jarowski, C.I. Effect of selected amino acids on ethanol toxicity in rats. *J. Pharm. Sci.* 62(1), 49-55 (1973).

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, 12/12/2022

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