

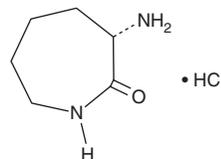
PRODUCT INFORMATION



L-Lysine lactam (hydrochloride)

Item No. 31685

CAS Registry No.: 26081-07-2
Formal Name: (3S)-3-aminohexahydro-2H-azepin-2-one, monohydrochloride
Synonyms: L-(-)- α -Amino- ϵ -caprolactam, (S)-3-Amino-2-azepanone
MF: C₆H₁₂N₂O • HCl
FW: 164.6
Purity: \geq 95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

L-lysine lactam (hydrochloride) is supplied as a crystalline solid. Aqueous solutions of L-lysine lactam (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of L-lysine 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-lysine lactam is a building block.^{1,2} It has been used in the synthesis of lysine sulfonamide HIV protease inhibitors, as well as bengamide derivatives with *in vitro* anticancer activity.

References

1. Stranix, B.R., Lavallée, J.-F., Sévigny, G., *et al.* Lysine sulfonamides as novel HIV-protease inhibitors: *N* ϵ -Acyl aromatic α -amino acids. *Bioorg. Med. Chem. Lett.* **16(13)**, 3459-3462 (2006).
2. Phi, T.D., Mai, H.D.T., Tran, V.H., *et al.* Design, synthesis and cytotoxicity of bengamide analogues and their epimers. *Medchemcomm.* **8(2)**, 445-451 (2016).

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, 13/09/2022

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