

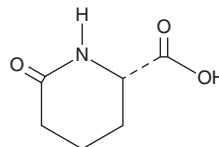
PRODUCT INFORMATION



L-Pyrohomo-glutamic Acid

Item No. 31694

CAS Registry No.: 34622-39-4
Formal Name: (2S)-6-oxo-2-piperidinecarboxylic acid
Synonyms: L-Pyro- α -aminoadipic Acid,
6-Oxo-L-pipecolic Acid
MF: C₆H₉NO₃
FW: 143.1
Purity: $\geq 95\%$
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

L-Pyrohomo-glutamic acid is supplied as a solid. A stock solution may be made by dissolving the L-pyrohomo-glutamic acid in the solvent of choice, which should be purged with an inert gas. L-Pyrohomo-glutamic acid is soluble in organic solvents such as ethanol and DMSO. The solubility of L-pyrohomo-glutamic acid in these solvents is approximately 0.2 and 1 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of L-pyrohomo-glutamic acid can be prepared by directly dissolving the L-pyrohomo-glutamic acid in aqueous buffers. The solubility of L-pyrohomo-glutamic acid 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-Pyrohomo-glutamic acid is an amino acid building block.^{1,2} It has been used in the synthesis of ligands for FK506-binding proteins (FKBPs) and histone deacetylase (HDAC) inhibitors.

References

1. Pomplun, S., Wang, Y., Kirschner, A., *et al.* Rational design and asymmetric synthesis of potent and neurotrophic ligands for FK506-binding proteins (FKBPs). *Angew. Chem. Int. Ed.* **54**(1), 345-348 (2015).
2. Taddei, M., Cini, E., Giannotti, L., *et al.* Lactam based 7-amino suberoylamide hydroxamic acids as potent HDAC inhibitors. *Bioorg. Med. Chem. Lett.* **24**(1), 61-64 (2014).

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, 12/09/2022

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