

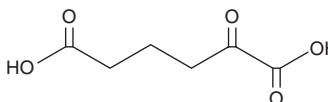
PRODUCT INFORMATION



2-Oxoadipic Acid

Item No. 34574

CAS Registry No.: 3184-35-8
Formal Name: 2-oxo-hexanedioic acid
Synonyms: α -Ketoadipate, 2-Ketoadipic Acid, α -Ketoadipic Acid, α -Oxoadipic Acid
MF: C₆H₈O₅
FW: 160.1
Purity: \geq 95%
Supplied as: A 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

2-Oxoadipic acid is supplied as a solid. A stock solution may be made by dissolving the 2-oxoadipic acid in water. The solubility of 2-oxoadipic acid in water (sonicated) is approximately 250 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

2-Oxoadipic acid is an intermediate in the catabolism of L-tryptophan, L-lysine, and hydroxylysine.^{1,2} Urinary excretion of 2-oxoadipic acid is increased in patients with α -ketoadipic aciduria, a rare inborn error in the metabolism of 2-oxoadipic acid to glutaryl-coenzyme A (glutaryl-CoA).

References

1. Xia, Z.-W., Inoue, Y., Ohse, M., *et al.* A study on α -ketoadipic aciduria by gas chromatographic-mass spectrometry. *World J. Gastroenterol.* **6(5)**, 766-769 (2000).
2. Gray, R.G., O'Neill, E.M., and Pollitt, R.J. α -aminoadipic aciduria: Chemical and enzymatic studies. *J. Inherit. Metab. Dis.* **2(4)**, 89-92 (1980).

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, 01/03/2023

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