

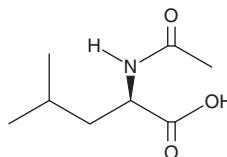
PRODUCT INFORMATION



N-acetyl-D-Leucine

Item No. 30489

CAS Registry No.: 19764-30-8
MF: $C_8H_{15}NO_3$
FW: 173.2
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

N-acetyl-D-Leucine is supplied as a crystalline solid. A stock solution may be made by dissolving the N-acetyl-D-leucine in the solvent of choice, which should be purged with an inert gas. N-acetyl-D-Leucine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of N-acetyl-D-leucine in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

N-acetyl-D-Leucine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, N-acetyl-D-leucine should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. N-acetyl-D-Leucine has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

N-acetyl-D-Leucine is a derivative of D-leucine and a substrate for various enzymes in the amidohydrolase superfamily.¹ It has been used to characterize the function of amidohydrolase enzymes from *B. bronchiseptica*, *S. coelicolor*, and *G. oxydans*.

Reference

1. Cummings, J., Fedorov, A.A., Xu, C., *et al.* Annotating enzymes of uncertain function: The deacylation of D-amino acids by members of the amidohydrolase superfamily. *Biochemistry* **48**(27), 6469-6481 (2009).

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, 12/01/2022

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