

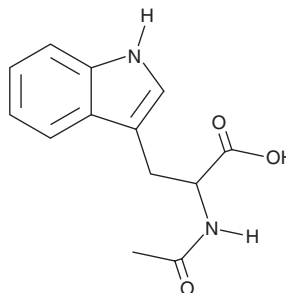
PRODUCT INFORMATION



N-acetyl-DL-Tryptophan

Item No. 36399

CAS Registry No.: 87-32-1
Formal Name: N-acetyl-tryptophan
Synonyms: DL-N-Acetyltryptophan, DL-NAT, NSC 49124
MF: C₁₃H₁₄N₂O₃
FW: 246.3
Purity: ≥98%
UV/Vis.: λ_{max}: 222, 282, 291 nm
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

N-acetyl-DL-Tryptophan (DL-NAT) is supplied as a solid. A stock solution may be made by dissolving the DL-NAT in the solvent of choice, which should be purged with an inert gas. DL-NAT is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of DL-NAT in these solvents is approximately 15 and 20 mg/ml, respectively. DL-NAT is slightly soluble in ethanol.

Description

DL-NAT is a derivative of the amino acid DL-tryptophan (Item No. 31210) and a racemic mixture of L-NAT and D-NAT.¹ It protects against hydrogen peroxide-induced cell death of NSC-34 motoneurons in an *in vitro* model of amyotrophic lateral sclerosis (ALS; IC₅₀ = 1 μM).

Reference

1. Sirianni, A.C., Jiang, J., Zeng, J., *et al.* N-acetyl-L-tryptophan, but not N-acetyl-D-tryptophan, rescues neuronal cell death in models of amyotrophic lateral sclerosis. *J. Neurochem.* **135**(5), 956-968 (2015).

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, 02/03/2023

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