

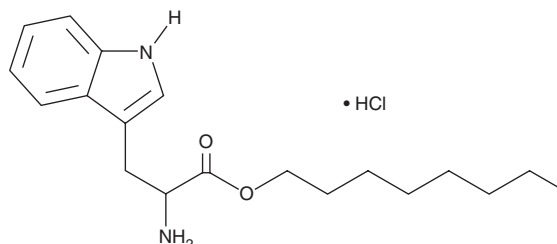
PRODUCT INFORMATION



DL-Tryptophan octyl ester (hydrochloride)

Item No. 17599

CAS Registry No.: 6278-90-6
Formal Name: tryptophan, octyl ester, monohydrochloride
Synonym: NSC 34501
MF: C₁₉H₂₈N₂O₂ • HCl
FW: 352.9
Purity: ≥98%
UV/Vis.: λ_{max}: 220, 280, 288 nm
Supplied as: A crystalline 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

DL-Tryptophan octyl ester (hydrochloride) is supplied as a crystalline solid, which should be purged with an inert gas. A stock solution may be made by dissolving the DL-tryptophan octyl ester (hydrochloride) in the solvent of choice. DL-Tryptophan octyl ester (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of DL-tryptophan octyl ester (hydrochloride) in these solvents is approximately 1, 25, and 15 mg/ml, respectively.

DL-Tryptophan octyl ester (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, DL-tryptophan octyl ester (hydrochloride) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. DL-Tryptophan octyl ester (hydrochloride) has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Tryptophan is an amino acid precursor of serotonin and melatonin. DL-Tryptophan octyl ester is a tryptophan derivative with an octyl ester protected COOH group.¹ This addition creates a stable, cell-permeable molecule capable of generating free tryptophan upon hydrolysis of the ester bond.

Reference

1. Nakamura, H. and Pisano, J.J. Sensitive and specific detection of tryptophan, tryptamine and N-terminal tryptophan peptides on thin-layer plates using a unique fluorogenic reaction with fluorescamine. *J. Chromatogr.* **152**, 153-165 (1978).

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, 11/10/2022

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