

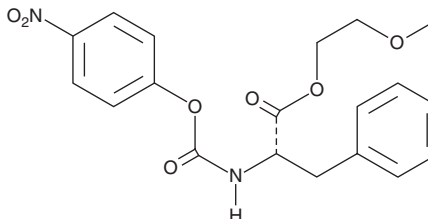
PRODUCT INFORMATION



(S)-NIFE

Item No. 17431

CAS Registry No.: 328406-65-1
Formal Name: N-[(4-nitrophenoxy)carbonyl]-L-phenylalanine, 2-methoxyethyl ester
MF: C₁₉H₂₀N₂O₇
FW: 388.4
Purity: ≥95%
UV/Vis.: λ_{max}: 275 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

(S)-NIFE is supplied as a crystalline solid. A stock solution may be made by dissolving the (S)-NIFE in the solvent of choice, which should be purged with an inert gas. (S)-NIFE is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of (S)-NIFE in these solvents is approximately 3, 10, and 20 mg/ml, respectively.

(S)-NIFE is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (S)-NIFE should first be dissolved in DMF and then diluted with the aqueous buffer of choice. (S)-NIFE has a solubility of approximately 0.12 mg/ml in a 1:7 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

(S)-NIFE is a chiral derivatizing agent for amino acid analysis by HPLC.¹ It can also be used for derivatization of secondary amino acids.¹

Reference

1. Péter, A., Vékes, E., Géza, T., *et al.* Application of a new chiral derivatizing agent to the enantioseparation of secondary amino acids. *J. Chromatogr. A.* **948(1-2)**, 283-294 (2002).

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, 12/08/2022

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