

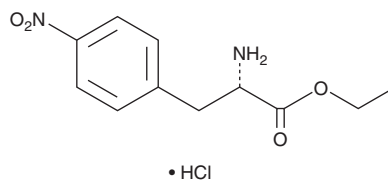
PRODUCT INFORMATION



Ethyl 4-nitro-L-phenylalanine (hydrochloride)

Item No. 30615

CAS Registry No.: 58816-66-3
Formal Name: 4-nitro-L-phenylalanine, ethyl ester, monohydrochloride
MF: C₁₁H₁₄N₂O₄ • HCl
FW: 274.7
Purity: ≥98%
UV/Vis.: λ_{max}: 266 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

Ethyl 4-nitro-L-phenylalanine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ethyl 4-nitro-L-phenylalanine should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Ethyl 4-nitro-L-phenylalanine has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Ethyl 4-nitro-L-phenylalanine is a building block.^{1,2} It has been used in the synthesis of chemokine (C-C motif) receptor 3 (CCR3) antagonists, as well as α4β1 (VLA-4) integrin antagonists.

References

1. White, J.R., Lee, J.M., Dede, K., *et al.* Identification of potent, selective non-peptide CC chemokine receptor-3 antagonist that inhibits eotaxin-, eotaxin-2-, and monocyte chemoattractant protein-1-induced eosinophil migration. *J. Biol. Chem.* **275**(47), 36626-36631 (2000).
2. Phillips, D.J., Davenport, R.J., Demaude, T.A., *et al.* Imidazopyridines as VLA-4 integrin antagonists. *Bioorg. Med. Chem. Lett.* **18**(14), 4146-4149 (2008).

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