

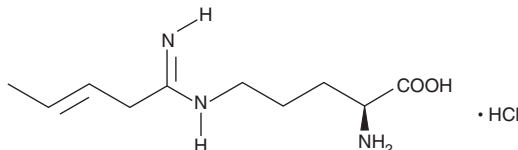
PRODUCT INFORMATION



Propenyl-L-NIO (hydrochloride)

Item No. 10011724

Formal Name: N⁵-[(3E)-1-imino-3-pentenyl]-L-ornithine, monohydrochloride
Synonym: ENIPO
MF: C₁₀H₁₉N₃O₂ • HCl
FW: 249.7
Purity: ≥90%
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

ENIPO (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the ENIPO (hydrochloride) in the organic solvent of choice, which should be purged with an inert gas. ENIPO (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of ENIPO (hydrochloride) in these solvents is approximately 10, 2, and 2.5 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of ENIPO (hydrochloride) can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of ENIPO (hydrochloride) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

ENIPO (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the ENIPO (hydrochloride) in an organic solvent purged with an inert gas. ENIPO (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ENIPO (hydrochloride) in ethanol is approximately 10 mg/ml and approximately 2 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of ENIPO (hydrochloride) can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of ENIPO (hydrochloride) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Reference

1. Bretscher, L.E., Li, H., Poulos, T.L., *et al.* Structural characterization and kinetics of nitric-oxide synthase inhibition by novel N⁵-(iminoalkyl)- and N⁵-(iminoalkenyl)-ornithines. *J. Biol. Chem.* **278**(47), 46789-46797 (2003).

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, 12/12/2022

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