

PRODUCT INFORMATION



NOC-5

Item No. 16534

CAS Registry No.: 146724-82-5

Formal Name: 3-[2-hydroxy-1-(1-methylethyl)-2-nitrosohydrazinyl]-1-propanamine

MF: C₆H₁₆N₄O₂

FW: 176.2

Purity: ≥98%

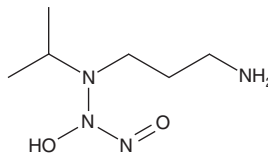
UV/Vis.: λ_{max}: 251 nm

Supplied as: A crystalline solid

Storage: -80°C

Stability: As supplied, 2 years from the QC date provided on the Certificate of Analysis, when stored properly

Special Conditions: Light, oxygen, and temperature sensitive



Laboratory Procedures

NOC-5 is supplied as a crystalline solid. NOC-5 is sparingly soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. For biological experiments, we suggest that organic solvent-free aqueous solutions of NOC-5 be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of NOC-5 in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

NOC-5 is a stable nitric oxide (NO)-amine complex that acts as a NO donor.¹ It can release two equivalents of NO in solution under physiological conditions without a cofactor.¹ The half-life of NOC-5 in PBS (pH 7.4) is 93 minutes at 22°C, and it remains relatively stable in alkaline solution (pH ≥10.0).¹

Reference

1. Hrabie, J.A., Klose, J.R., Wink, D.A., *et al.* New nitric oxide-releasing zwitterions derived from polyamines. *J. Org. Chem.* **58**, 1472-1476 (1993).

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, 05/23/2016

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