

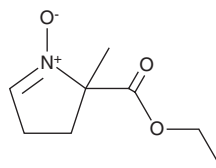
PRODUCT INFORMATION



EMPO

Item No. 20618

CAS Registry No.: 61856-99-3
Formal Name: 3,4-dihydro-2-methyl-2H-Pyrrole-2-carboxylic acid 1-oxide, ethyl ester
Synonym: 5-(Ethoxycarbonyl)-5-methyl-1-Pyrroline-N-Oxide
MF: C₈H₁₃NO₃
FW: 171.2
Purity: ≥95%
UV/Vis.: λ_{max}: 238 nm
Supplied as: A 25 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

EMPO is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of EMPO in these solvents is approximately 15 and 30 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. If an organic solvent-free solution of EMPO is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of EMPO in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

EMPO is a hydrophilic cyclic nitron analog of the free radical spin trap DMPO (Item No. 10006436). The EMPO-superoxide adduct (EMPO-OOH) exhibits an ESR spectrum that is similar to the DMPO-OOH spectrum.¹ However, the EMPO-OOH adduct is about 5-8 times more stable ($t_{1/2}$ = 4.8 – 8.6 min) and does not spontaneously decay to the hydroxyl adduct, making spectral interpretation less difficult.¹⁻³

References

1. Olive, G., Mercier, A., Le Moigne, F., *et al.* 2-ethoxycarbonyl-2-methyl-3,4-dihydro-2H-pyrrole-1-oxide: Evaluation of the spin trapping properties. *Free Radic. Biol. Med.* **28**(3), 403-408 (2000).
2. Stolze, K., Udilova, N., and Nohl, H. Spin adducts of superoxide, alkoxyl, and lipid-derived radicals with EMPO and its derivatives. *Biol. Chem.* **383**(5), 813-820 (2002).
3. Zhang, H., Joseph, J., Vasquez-Vivar, J., *et al.* Detection of superoxide anion using an isotopically labeled nitron spin trap: Potential biological applications. *FEBS Lett.* **473**(1), 58-62 (2000).

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, 12/18/2025

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