

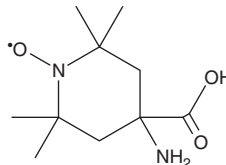
PRODUCT INFORMATION



TOAC

Item No. 42097

CAS Registry No.: 15871-57-5
Formal Name: 4-amino-4-carboxy-2,2,6,6-tetramethyl-1-piperidinyloxy
Synonym: NSC 239127
MF: C₁₀H₁₉N₂O₃
FW: 215.3
Purity: ≥95%
Supplied as: A 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

TOAC is supplied as a solid. Aqueous solutions of TOAC can be prepared by directly dissolving the solid in aqueous buffers. TOAC is slightly soluble (0.1-1 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

TOAC is a nitroxyl amino acid and spin label.^{1,2} It incorporates a stable nitrogen radical into peptide backbones, which allows it to function as an intrinsic reporter during electron paramagnetic resonance (EPR) spectroscopy, and has been used to study peptide structural dynamics.

References

1. Corvaja, C., Sartori, E., Toffoletti, A., *et al.* Interaction between TOAC free radical and photoexcited triplet chromophores linked to peptide templates. *Biopolymers* **55(6)**, 486-495 (2000).
2. Marsh, D. Orientation of TOAC amino-acid spin labels in α -helices and β -strands. *J. Magn. Reson.* **180(2)**, 305-310 (2006).

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, 02/04/2025

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