

Carboxy-PTIO

Nitric oxide (NO) scavenger

Carboxy-PTIO is a water-soluble and stable nitric oxide radical scavenger that shows in chemical and biological systems antagonistic action against the free nitric oxide radical ($\text{NO}\cdot$) in a stoichiometric manner. The inhibitory effects are almost two-fold stronger than those of NMMA (Prod. No. ALX-106-001) and NOARG (Prod. No. ALX-105-001). Carboxy-PTIO can be used for an improved detection of nitric oxide in activated macrophage cultures in combination with the Griess reagent (Prod. No. ALX-400-004). Carboxy-PTIO also protects against endotoxin shock.

Citations: 17

[View Online »](#)

Ordering Information

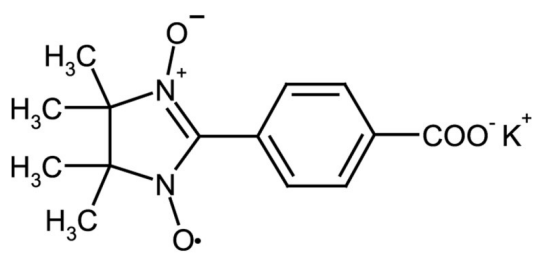
[Order Online »](#)

ALX-430-001-M010	10mg
ALX-430-001-M050	50mg
ALX-430-001-M250	250mg

Manuals, SDS & CofA

[View Online »](#)

- Highly pure NO scavenger



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 2-(4-Carboxyphenyl)-4,4,5,5-tetramethylimidazoline-1-oxyl-3-oxide . potassium salt

Appearance Blue solid.

CAS 148819-94-7

Couple Target Nitric oxide

Formula $C_{14}H_{16}N_2O_4 \cdot K$

Identity Identity determined by MS and NMR.

MW 315.4

Purity ≥99% (HPLC)

Solubility Soluble in DMF, DMSO, 100% ethanol, or 100% methanol. Highly soluble in water or aqueous buffers (e.g. PBS, pH 7.2).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)