# **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 (NO·) 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**

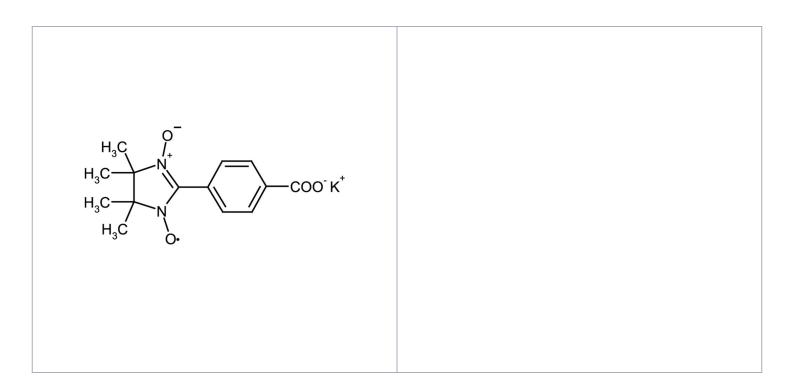
**Order Online** »

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  $C_{14}H_{16}N_2O_4$  . 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).

Last modified: May 29, 2024

