

# GEA 5024

## Nitric oxide donor

Water soluble and stable nitric oxide donor.

Citations: 4

[View Online »](#)

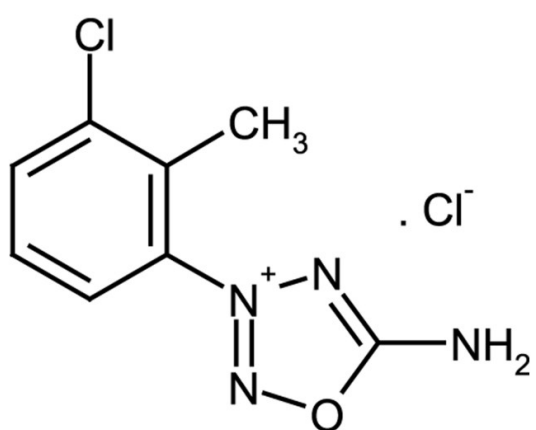
### Ordering Information

[Order Online »](#)

|                  |      |
|------------------|------|
| ALX-430-005-M010 | 10mg |
|------------------|------|

### Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

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

**Long Term Storage** +4°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 5-Amino-3-(3-chloro-2-methylphenyl)-1,2,3,4-oxatriazolium chloride

**Appearance** White to off-white crystalline solid.

**CAS** 144575-27-9

**Formula**  $C_8H_7ClN_4O \cdot HCl$

**MW** 210.6 . 36.5

**Purity** ≥97% (HPLC)

**Solubility** Soluble in water.

**Last modified: May 29, 2024**



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-  
be@enzolifesciences.com](mailto:info-be@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)