

# GEA 3162

## Nitric oxide donor

Water soluble nitric oxide donor and potent inhibitor of ADP-induced platelet aggregation in platelet rich plasma (PRP). Stimulates cGMP production in platelets, granulocytes and polymorphonuclear leukocytes. Vasodepressant.

Citations: 11

[View Online »](#)

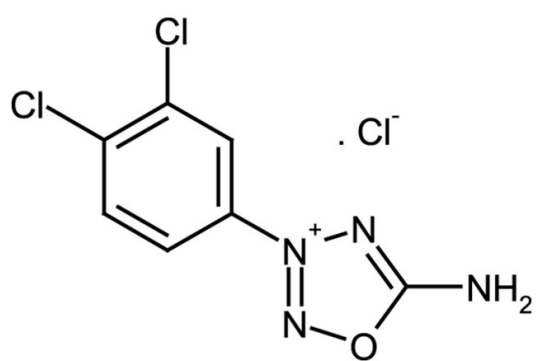
## Ordering Information

[Order Online »](#)

|                  |      |
|------------------|------|
| ALX-430-004-M010 | 10mg |
|------------------|------|

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

|                   |                                                                  |
|-------------------|------------------------------------------------------------------|
| Use/Stability     | As indicated on product label or CoA when stored as recommended. |
| Handling          | Protect from oxygen. Use fresh solutions.                        |
| Long Term Storage | -20°C                                                            |
| Shipping          | Ambient Temperature                                              |

## Regulatory Status

RUO - Research Use Only

## Product Details

|                  |                                                                  |
|------------------|------------------------------------------------------------------|
| Alternative Name | 1,2,3,4-Oxatriazolium, 5-amino-3-(3,4-dichlorophenyl)-, chloride |
| Appearance       | White to off-white crystalline solid.                            |
| CAS              | 144575-47-3                                                      |
| Formula          | $C_7H_4Cl_2N_4O \cdot HCl$                                       |
| Identity         | Determined by NMR.                                               |
| MW               | 231.0 . 36.5                                                     |
| Purity           | ≥97% (TLC)                                                       |
| Solubility       | Soluble in water, methanol or DMSO.                              |

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-  
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)