

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 »](#)

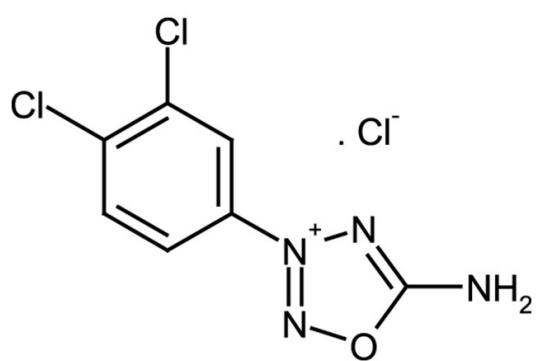
Ordering Information

[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

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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