

# NS-2028

## Guanylyl cyclase inhibitor

A potent and selective, cell-permeable inhibitor of nitric oxide (NO)-sensitive soluble guanylyl cyclase (sGC). Inhibition is irreversible. Inhibits also the basal, the YC-1 and the coenhanced sGC activity. Complete inhibition of NO-induced relaxation in vascular and nonvascular smooth muscle at 1µM.

Citations: 5

[View Online »](#)

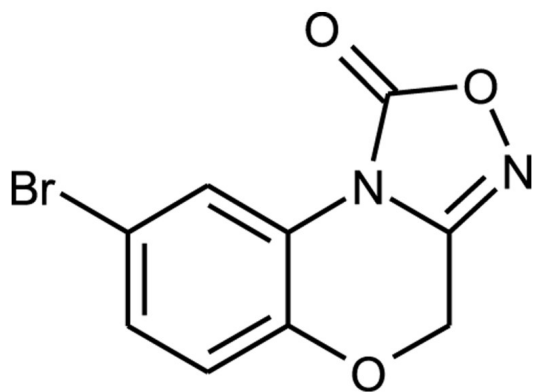
## Ordering Information

[Order Online »](#)

ALX-270-210-M001	1mg
ALX-270-210-M005	5mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended.
<b>Handling</b>	Protect from moisture.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	4H-8-Bromo-1,2,4-oxadiazolo(3,4-d)benz(b)(1,4)oxazin-1-one
<b>Appearance</b>	White to off-white crystalline powder.
<b>CAS</b>	204326-43-2
<b>Couple Target</b>	Guanylyl cyclase
<b>Couple Type</b>	Inhibitor
<b>Formula</b>	$C_9H_5BrN_2O_3$
<b>MW</b>	269.1
<b>Purity</b>	≥98% (TLC)
<b>Solubility</b>	Soluble in DMSO or 100% ethanol; insoluble 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)