

# LY-83,583

## Guanylyl cyclase inhibitor

Competitive inhibitor of soluble guanylyl cyclase (sGC) ( $IC_{50}=2\mu M$ ) and of cGMP production. Neurotoxin. Blocks antigen-induced leukotriene release and reduces platelet-endothelial cell adhesion. Blocks acetylcholine-induced vasorelaxation.

Citations: 12

[View Online »](#)

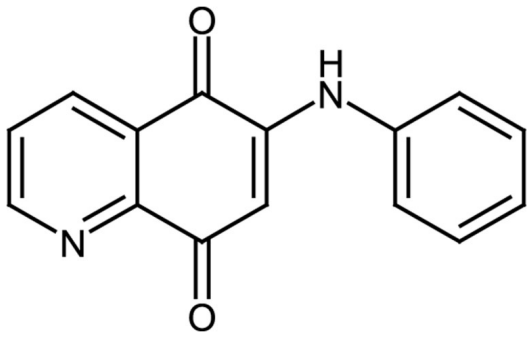
## Ordering Information

[Order Online »](#)

ALX-550-002-M005	5mg
ALX-550-002-M025	25mg

## Manuals, SDS & CofA

[View Online »](#)



## 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** 6-(Phenylamino)-5,8-quinolinedione, 6-Anilino-5,8-quinolinedione

**Appearance** Violet to dark red powder.

**CAS** 91300-60-6

**Couple Target** Guanylyl cyclase

**Couple Type** Inhibitor

**Formula**  $C_{15}H_{10}N_2O_2$

**Identity** Determined by <sup>1</sup>H-NMR.

**MW** 250.3

**Purity** ≥95% (TLC)

**Solubility** Soluble in methanol, methylene chloride or diluted aqueous acid; insoluble in water.

**Source** Synthetic.

**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)