

# Furoxan

Releases NO in the presence of thiols

A novel NO donor which only releases NO in the presence of thiols. It is a potent vasodilator and inhibitor of platelet aggregation.

Citations: 2

[View Online »](#)

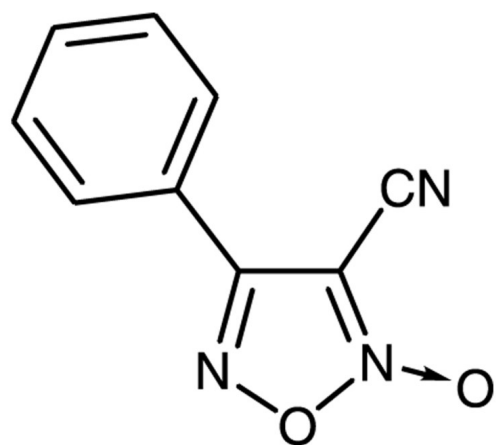
Ordering Information

[Order Online »](#)

BML-CN255-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	4-Phenyl-3-furoxanarbonitrile
<b>Appearance</b>	White solid.
<b>CAS</b>	125520-62-9
<b>Formula</b>	$C_9H_5N_3O_2$
<b>MW</b>	187.2
<b>MeltingPoint</b>	72-74°C
<b>Purity</b>	≥99% (HPLC)
<b>Solubility</b>	Soluble in ethanol (25 mg/ml) or DMSO (25 mg/ml).

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)