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 »

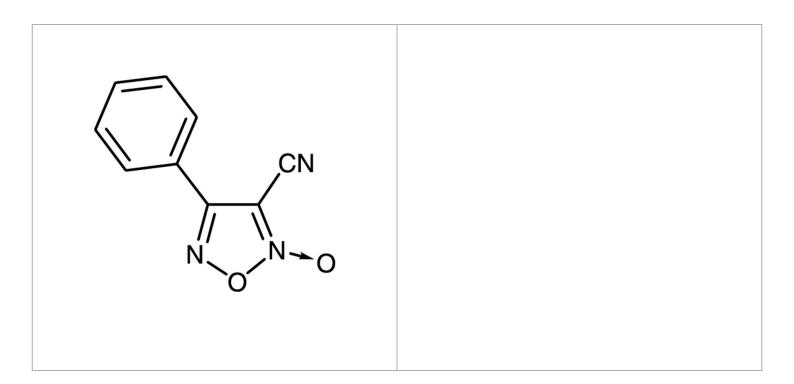
Ordering Information

Order Online »

BML-CN255-0010	10mg
BML-CN255-0050	50mg

Manuals, SDS & CofA

View Online »



Handling & Storage

Use/Stability 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.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 4-Phenyl-3-furoxancarbonitrile

Appearance White solid.

CAS 125520-62-9

Formula $C_9H_5N_3O_2$

MW 187.2

MeltingPoint 72-74°C

Purity ≥99% (HPLC)

Solubility Soluble in ethanol (25 mg/ml) or DMSO (25 mg/ml).

Last modified: May 29, 2024



eu@enzolifesciences.com