

# SNAP

## Releases NO under physiological conditions

Stable analog of endogenous S-nitroso compounds. Nitric oxide (NO) donor and a source of NO *in vivo* which spontaneously releases NO under physiological conditions. The half-life time is 4.6 hours in aqueous media. Potent vasodilator *in vitro* and *in vivo* that does not induce pharmacological tolerance. Relaxes isolated bovine coronary artery rings with an  $EC_{50}=0.1\mu M$ . It also inhibits vascular smooth muscle cell mitogenesis and proliferation, inhibits leukocyte adherence to endothelium, and inhibits cysteine proteases. The unique NO releasing properties of SNAP make it an ideal agent for studying the pharmacological and physiological actions of NO.

Citations: 20

[View Online »](#)

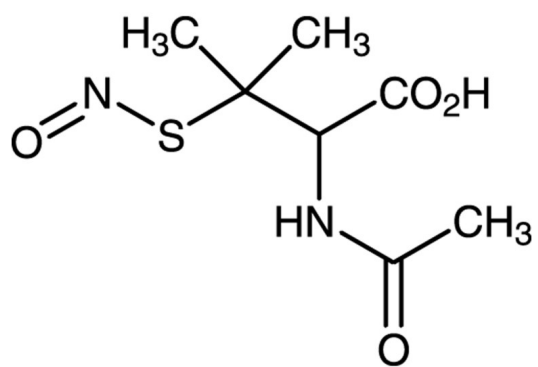
## Ordering Information

[Order Online »](#)

BML-CN210-0020	20mg
BML-CN210-0100	100mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for up to 1 year after receipt when stored at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	S-Nitroso-N-acetyl-D,L-penicillamine
Appearance	Green crystalline solid.
CAS	79032-48-7
Couple Target	Protease
Couple Type	Inhibitor
Formula	$C_7H_{12}N_2O_4S$
MW	220.3
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (25mg/ml) 100% ethanol (25mg/ml) or methanol; slightly soluble in water (2mg/ml).
Source	Synthetic.
Technical Info / Product Notes	Note: Product is not sterile.

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-fr@enzolifesciences.com](mailto:info-fr@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)