

# 2-Iminobiotin

NOS inhibitor

Reversible inhibitor of mouse inducible and rat neuronal nitric oxide synthase (iNOS/NOS II; nNOS/NOS I).

Citations: 1

[View Online »](#)

## Ordering Information

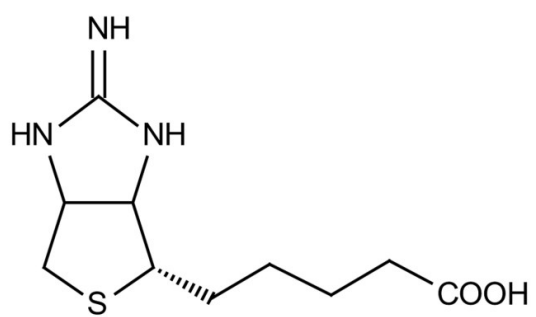
[Order Online »](#)

ALX-430-008-M010

10mg

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** +4°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Guanidinobiotin

**Appearance** White to off-white powder.

**CAS** 13395-35-2

**Couple Target** NOS

**Couple Type** Inhibitor

**Formula**  $C_{10}H_{17}N_3O_2S$

**Identity** Determined by IR.

**MW** 243.3

**Purity** ≥98% (TLC)

**Solubility** Soluble in 0.1N hydrochloric acid.

**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)