

SIN-1A/ γ CD Complex

Nitric oxide donor

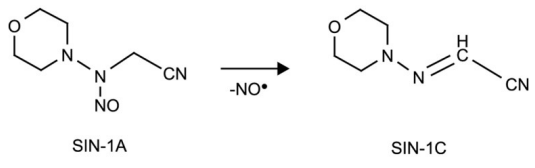
Physiologically active nitric oxide (NO) releasing agent. Releasing one equivalent of NO the half life time at room temperature is ~40 minutes mostly independent from the pH.

Citations: 2

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	3-Morpholino-N-nitroso-aminoacetonitrile . γ -cyclodextrin complex
Appearance	Pale yellow powder.
CAS	26687-79-6 (SIN-1A)
Formulation	The complex contains $10 \pm 2\%$ of SIN-1A. SIN-1 content: not detectable. SIN-1C (N-morpholinoiminoacetonitrile) content: $0.5 \pm 0.2\%$.
Quantity	* Sold by net SIN-1A content.
Solubility	Well soluble in distilled water, isotonic saline, phosphate buffer or HCl solution, pH 1.4.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com