

Guanylyl cyclase (soluble) (bovine lung)

Citations: 4

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Product should be active for at least one year after shipping date when stored at -80°C under nitrogen in the presence of 50% glycerol. Dilutions of the enzyme should be performed in the presence of 0.5mg/ml BSA. The dissolved product is stable for 8 months when stored in 50% glycerol at -80°C.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	GTP pyrophosphate-lyase (cycling)
Concentration	
Formulation	Liquid. In ~20µl containing 50% glycerol, 88mM NaCl, 1mM DTT, 0.5mM EDTA, 25mM triethanolamine at pH 7.4.
Source	Isolated from bovine lung by an immunoaffinity method.
Specific Activity	>10'000nmol cGMP/mg/min at 37°C under NO-stimulated conditions.

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)