

Guanylin (rat, mouse)

Guanylyl cyclase activator

Endogenous intestinal guanylyl cyclase activator.

Intestinal guanylate cyclase possesses an extracellular ligand binding domain, a transmembrane domain, an intracellular protein kinase-like region and a cyclase catalytic domain. GC-C is expressed in various mammalian tissues including intestine, adrenal gland, airway epithelial cells, brain, and olfactory and tracheal mucosa.

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-157-011-MC01	0.1mg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance White to off-white powder.

CAS 144940-98-7

Formula $C_{60}H_{90}N_{16}O_{22}S_4$

Formulation Lyophilized

MW 1515.7

Purity ≥99%

Sequence H-Pro-Asn-Thr-Cys-Glu-Ile-Cys-Ala-Tyr-Ala-Ala-Cys-Thr-Gly-Cys-OH(Disulfide bonds between Cys⁴-Cys¹² and Cys⁷-Cys¹⁵).

Solubility Soluble in water.

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