

Protoporphyrin IX (free acid)

Guanylyl cyclase activator

Protoporphyrin IX free acid, as distinct from its zinc salt (which inhibits heme oxygenase), activates soluble guanylyl cyclase (sGC) by binding directly to the enzyme. A useful reagent in cases where the use of nitric oxide (NO) or nitric oxide donors is undesirable.

Citations: 8

[View Online »](#)

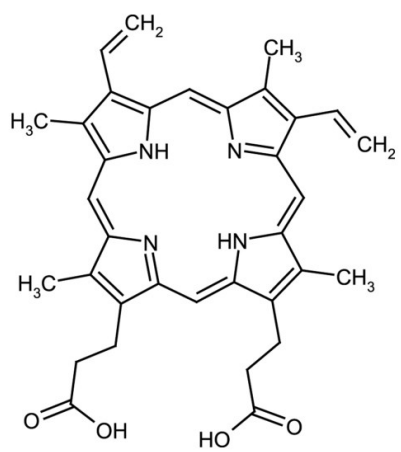
Ordering Information

[Order Online »](#)

ALX-430-041-M300	300mg
ALX-430-041-G001	1g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Make solutions fresh daily.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	8,13-Divinyl-3,7,12,17-tetramethyl-21H,23H-porphine-2,18-dipropionic acid
Appearance	Purple to dark brown solid.
CAS	553-12-8
Couple Target	Guanylyl cyclase
Couple Type	Activator
Formula	$C_{34}H_{34}N_4O_4$
Identity	Determined by NMR and UV/VIS.
MI	14: 7897
MW	562.7
Purity	≥95% (HPLC)
Solubility	Dissolve initially in 0.1M base (Tris base or NaOH), then add water miscible organic solvent (EtOH (10mg/ml), MeOH, DMSO (10mg/ml), DMF) until solution is 50/50 base/solvent. When a clear solution is obtained, it can be diluted into an aqueous medium and titrated or buffered to any pH>7. Use immediately.



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)