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 »

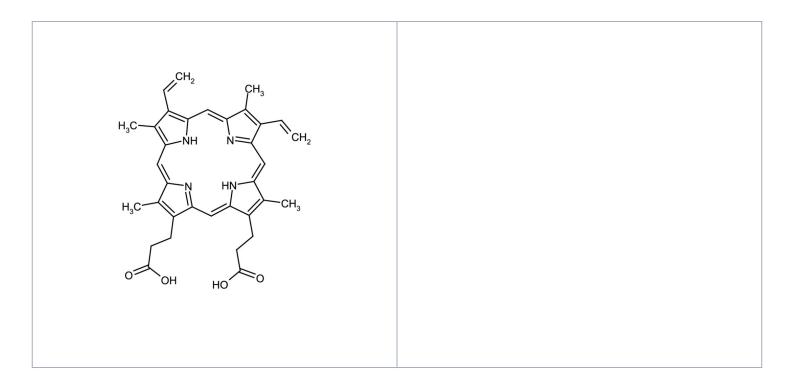
Ordering Information

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.

Last modified: May 29, 2024



info-

eu@enzolifesciences.com