

YM-254890

Gq inhibitor

Cyclic depsipeptide from *Chromobacterium* sp. YM-254890 is a novel Gq/11 inhibitor, U46619-induced platelet aggregation is inhibited but phorbol ester- or ionophore-induced aggregation is not. It blocks *Pasteurella multocida* toxin-mediated activation of Gq. A highly useful new tool for studying Gq/11-coupled receptor signaling and resulting cellular events.

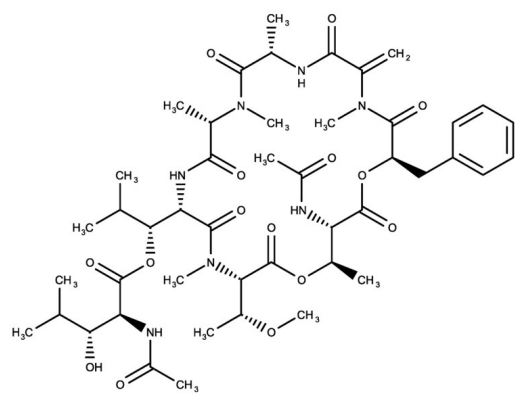
Ordering Information

[Order Online »](#)

ENZ-CHM486-0001	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 or off-white lyophilized solid.

CAS 568580-02-9

Couple Target G Protein

Couple Type Inhibitor

Formula $C_{46}H_{69}N_7O_{15}$

Identity Determined by NMR.

MW 960.10

Purity ≥95% (HPLC)

Solubility Soluble in DMSO (10 mg/ml).



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-
fr@enzolifesciences.com](mailto:info-fr@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)