

# 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.

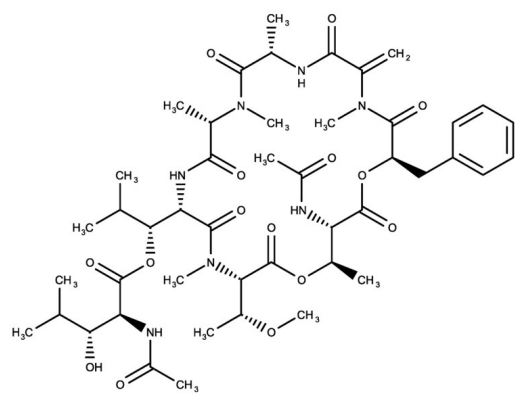
### Ordering Information

[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).

**Last modified: February 27, 2025**



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)