

# GP antagonist-2A

## G protein antagonist

A G protein antagonist which selectively inhibits the activation of  $G_q$  by  $M_1$  muscarinic cholinergic receptors (mAChR). Not effective in inhibiting  $M_2$  mAChR-promoted GTP hydrolysis by  $G_i$ . Inhibition is reversible and competitive with respect to receptor binding to G proteins. Also inhibits activation by cholecystokinin receptor in pancreatic acini at  $10\mu\text{M}$ .

Citations: 2

[View Online »](#)

## Ordering Information

[Order Online »](#)

BML-G515-0001	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** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** G-protein antagonist 2A

**Appearance** White to off-white solid.

**Formula**  $C_{79}H_{105}N_{21}O_{13}S$

**Formulation** Lyophilized.

**MW** 1588.9

**Purity** ≥99% (HPLC)

**Sequence** H-Arg-Pro-Lys-Pro-Gln-Gln-D-Trp-Phe-D-Trp-D-Trp-Met-NH<sub>2</sub>

**Solubility** Soluble in water (25mg/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-  
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)