

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-NH₂

Solubility Soluble in water (25mg/ml).

Last modified: May 29, 2024



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)