

# GP antagonist-2

## G protein antagonist

Selective, competitive and reversible G-protein antagonist. Inhibits the activation of  $G_i$  or  $G_o$  by  $M_2$  muscarinic cholinergic receptors (mAChR) and  $G_s$  by  $\beta$ -adrenergic receptors in reconstituted phospholipid vesicles. Inhibition is reversible and competitive with respect to receptor binding to G proteins. Substance P (Prod. No. ALX-167-003) related compound.

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	G-protein antagonist 2
<b>Appearance</b>	White to off-white solid.
<b>Formula</b>	$C_{57}H_{64}N_{12}O_9S$
<b>Formulation</b>	Lyophilized.
<b>MW</b>	1093.3
<b>Purity</b>	≥95% (HPLC)
<b>Sequence</b>	H-Pyr-Gln-D-Trp-Phe-D-Trp-D-Trp-Met-NH <sub>2</sub>
<b>Solubility</b>	Soluble in water or DMSO.
<b>Source</b>	Synthetic.

**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)