

SCH-202676

Blocks agonist and antagonist binding to G protein-coupled receptors

Modulator of both agonist and antagonist binding to G protein-coupled receptors which functions via thiol modification rather than true allosteric mechanisms. Has been shown to block agonist or antagonist activity at adenosine, adrenergic, dopaminergic, muscarinic, opioid and P2Y. IC_{50} s vary from 0.1 to 1.8 μ M. The nature of the interaction of SCH-202676 and receptors is dependent on whether it is studied using intact versus broken cell preparations.

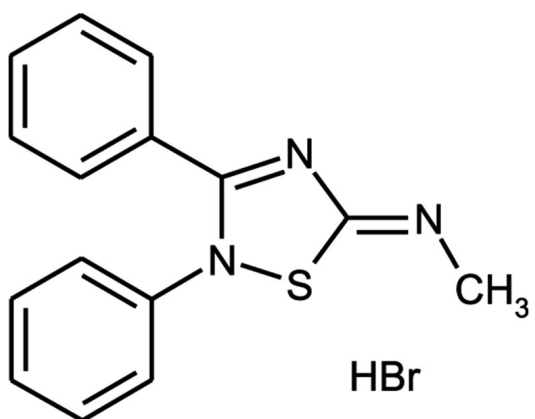
Ordering Information

[Order Online »](#)

BML-G521-0010	10mg
---------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance Yellow solid.

CAS 70375-43-8

Couple Target GPCR

Couple Type Blocker

Formula $C_{15}H_{13}N_3S \cdot HBr$

Identity Determined by NMR.

MW 348.3

Purity ≥99% (HPLC)

Solubility Soluble in DMSO (7mg/ml, with warming).



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)