# 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 $\mu$ M.The nature of the interaction of SCH-202676 and receptors is dependent on whether it is studied using intact versus broken cell preparations.

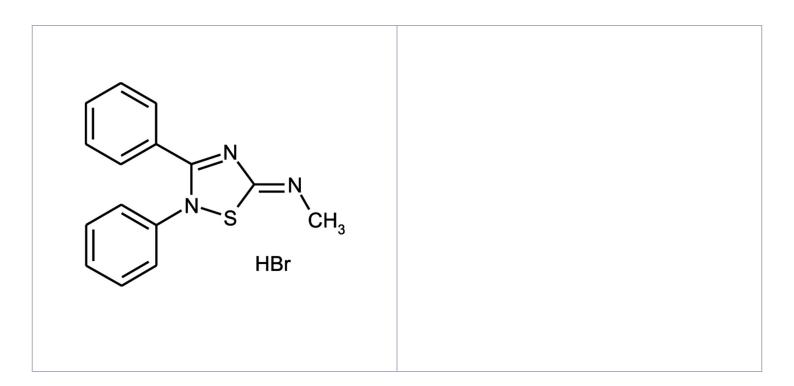
### **Ordering Information**

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  $C_{15}H_{13}N_3S$  . HBr

**Identity** Determined by NMR.

MW 348.3

Purity ≥99% (HPLC)

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

Last modified: May 29, 2024