

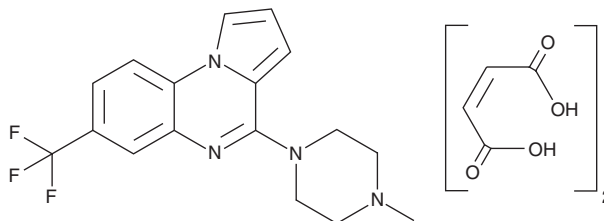
# PRODUCT INFORMATION



## CGS 12066B

Item No. 30623

**CAS Registry No.:** 109028-10-6  
**Formal Name:** 4-(4-methyl-1-piperazinyl)-7-(trifluoromethyl)-pyrrolo[1,2-a]quinoxaline, (2Z)-2-butenedioate  
**MF:** C<sub>17</sub>H<sub>17</sub>F<sub>3</sub>N<sub>4</sub> • 2C<sub>4</sub>H<sub>4</sub>O<sub>4</sub>  
**FW:** 566.5  
**Purity:** ≥98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

CGS 12066B is supplied as a solid. A stock solution may be made by dissolving the CGS 12066B in the solvent of choice, which should be purged with an inert gas. CGS 12066B is soluble in the organic solvent DMSO at a concentration of approximately 25 mg/ml.

### Description

CGS 12066B is an agonist of the serotonin (5-HT) receptor subtype 5-HT<sub>1B</sub> (IC<sub>50</sub> = 51 nM).<sup>1</sup> It is selective for 5-HT<sub>1B</sub> over 5-HT<sub>1A</sub> and 5-HT<sub>2</sub> receptors (IC<sub>50</sub>s = 876 and 6,480 nM, respectively), as well as α<sub>1</sub>-, α<sub>2</sub>-, and β-adrenergic, and dopamine D<sub>1</sub> and D<sub>2</sub> receptors (IC<sub>50</sub>s = >6,000, >1,000, >1,000, >5,000, and >5,000 nM, respectively). CGS 12066B inhibits forskolin-induced cAMP accumulation in opossum kidney cells when used at concentrations ranging from 0.001 to 1 μM.<sup>2</sup> *In vivo*, CGS 12066B reduces 5-hydroxy-L-tryptophan (5-HTP) accumulation in the fronto-parietal cortex (ED<sub>50</sub> = 7.92 μmol/kg) and induces dorsal raphe cell firing in rats (ED<sub>50</sub> = 358 nmol/kg).<sup>1</sup> It decreases interfemale and intermale aggression in Syrian hamsters.<sup>3</sup>

### References

1. Neale, R.F., Fallon, S.L., Boyar, W.C., *et al.* Biochemical and pharmacological characterization of CGS 12066B, a selective serotonin-1B agonist. *Eur. J. Pharmacol.* **136(1)**, 1-9 (1987).
2. Unsworth, C.D. and Molinoff, P.B. Regulation of the 5-hydroxytryptamine<sub>1B</sub> receptor in opossum kidney cells after exposure to agonists. *Mol. Pharmacol.* **42(3)**, 464-470 (1992).
3. Joppa, M.A., Rowe, R.K., and Meisel, R.L. Effects of serotonin 1A or 1B receptor agonists on social aggression in male and female Syrian hamsters. *Pharmacol. Biochem. Behav.* **58(2)**, 349-353 (1997).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/12/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM