

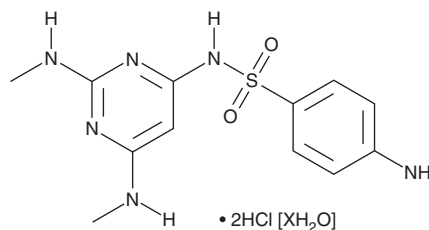
# PRODUCT INFORMATION



## Ro 04-6790 (hydrochloride hydrate)

Item No. 34476

**Formal Name:** 4-amino-N-[2,6-bis(methylamino)-4-pyrimidinyl]-benzenesulfonamide, dihydrochloride hydrate  
**MF:** C<sub>12</sub>H<sub>16</sub>N<sub>6</sub>O<sub>2</sub>S • 2HCl [XH<sub>2</sub>O]  
**FW:** 381.3  
**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

Ro 04-6790 (hydrochloride hydrate) is supplied as a solid. A stock solution may be made by dissolving the Ro 04-6790 (hydrochloride hydrate) in the solvent of choice, which should be purged with an inert gas. Ro 04-6790 (hydrochloride hydrate) is soluble in the organic solvent DMSO at a concentration of approximately 15 mg/ml. It is also soluble in water. The solubility of Ro 04-6790 (hydrochloride hydrate) in water is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Ro 04-6790 is an antagonist of the serotonin (5-HT) receptor subtype 5-HT<sub>6</sub> (K<sub>i</sub> = 44.7 nM).<sup>1</sup> It is selective for 5-HT<sub>6</sub> receptors over a panel of 23 additional receptors at 10 μM. Ro 04-6790 inhibits 5-HT-induced cAMP production in HeLa cells expressing human 5-HT<sub>6</sub> (pA<sub>2</sub> = 6.75). *In vivo*, Ro 04-6790 (10 and 30 mg/kg) induces stretching and yawning behavior in rats. It reverses scopolamine-induced decreases in novel object discrimination in rats.<sup>2</sup> Ro 04-6790 (3, 10, and 30 mg/kg) also attenuates cue-induced cocaine-seeking behavior in rats.<sup>3</sup>

### References

1. Sleight, A.J., Boess, F.G., Bös, M., *et al.* Characterization of Ro 04-6790 and Ro 63-0563: Potent and selective antagonists at human and rat 5-HT<sub>6</sub> receptors. *Br. J. Pharmacol.* **124**(3), 556-562 (1998).
2. Woolley, M.L., Marsden, C.A., Sleight, A.J., *et al.* Reversal of a cholinergic-induced deficit in a rodent model of recognition memory by the selective 5-HT<sub>6</sub> receptor antagonist, Ro 04-6790. *Psychopharmacology (Berl)* **170**(4), 358-367 (2003).
3. van Gaalen, M.M., Schettters, D., Schoffemeer, A.N.M., *et al.* 5-HT<sub>6</sub> antagonism attenuates cue-induced relapse to cocaine seeking without affecting cocaine reinforcement. *Int. J. Neuropsychopharmacol.* **13**(7), 961-965 (2010).

#### 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, 11/21/2022

#### 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