

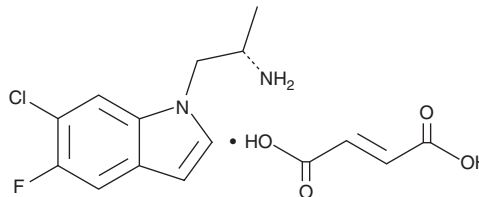
# PRODUCT INFORMATION



Ro 60-0175

Item No. 29520

**CAS Registry No.:** 169675-09-6  
**Formal Name:** (αS)-6-chloro-5-fluoro-α-methyl-1H-indole-1-ethanamine, (2E)-2-butenedioate  
**MF:** C<sub>11</sub>H<sub>12</sub>ClFN<sub>2</sub> • C<sub>4</sub>H<sub>4</sub>O<sub>4</sub>  
**FW:** 342.8  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 223, 280 nm  
**Supplied as:** A crystalline 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 60-0175 is supplied as a crystalline solid. A stock solution may be made by dissolving the Ro 60-0175 in the solvent of choice, which should be purged with an inert gas. Ro 60-0175 is soluble in the organic solvent DMSO.

## Description

Ro 60-0175 is an agonist of the serotonin (5-HT) receptor subtype 5-HT<sub>2</sub> that stimulates increases in intracellular calcium levels in CHO-K1 cells expressing human 5-HT<sub>2A</sub>, 5-HT<sub>2B</sub>, and 5-HT<sub>2C</sub> receptors (EC<sub>50</sub>s = 447, 0.9, and 32 nM, respectively).<sup>1</sup> It is selective for human 5-HT<sub>2A</sub> and 5-HT<sub>2C</sub> (K<sub>i</sub>s = 32 and 1 nM, respectively) over human 5-HT<sub>1A</sub>, 5-HT<sub>1D</sub>, 5-HT<sub>6</sub>, and 5-HT<sub>7</sub>, rat 5-HT<sub>1B</sub> and 5-HT<sub>3</sub>, and guinea pig 5-HT<sub>4</sub> receptors (K<sub>i</sub>s = ≥1,995 nM), as well as adenosine A<sub>1</sub> and A<sub>2</sub>, histamine H<sub>1-3</sub>, dopamine D<sub>1-5</sub>, α<sub>1</sub>-, α<sub>2</sub>-, and β<sub>1</sub>-adrenergic, M<sub>1-5</sub> muscarinic, nicotinic, kainate, AMPA, and μ-, δ-, and κ-opioid receptors (IC<sub>50</sub>s = >1,000 nM), but does bind to β<sub>2</sub>-adrenergic receptors (IC<sub>50</sub> = 251 nM).<sup>2</sup> *In vivo*, Ro 60-0175 increases plasma levels of corticosterone, oxytocin, and prolactin in rats (ED<sub>50</sub>s = 2.43, 4.19, and 4.03 mg/kg, respectively), effects that are not inhibited by the 5-HT<sub>2C</sub> antagonist SB-242084 (Item No. 10096) or the 5-HT<sub>2A</sub> antagonist MDL 100,907 (volinanserin; Item No. 15936).<sup>3</sup>

## References

- Porter, R.H., Benwell, K.R., Lamb, H., *et al.* Functional characterization of agonists at recombinant human 5-HT<sub>2A</sub>, 5-HT<sub>2B</sub> and 5-HT<sub>2C</sub> receptors in CHO-K1 cells. *Br. J. Pharmacol.* **128(1)**, 13-20 (1999).
- Martin, J.R., Bös, M., Jenck, F., *et al.* 5-HT<sub>2C</sub> receptor agonists: Pharmacological characteristics and therapeutic potential. *J. Pharmacol. Exp. Ther.* **286(2)**, 913-924 (1998).
- Damjanoska, K.J., Muma, N.A., Zhang, Y., *et al.* Neuroendocrine evidence that (S)-2-(chloro-5-fluoro-indol-1-yl)-1-methylethylamine fumarate (Ro 60-0175) is not a selective 5-hydroxytryptamine<sub>2C</sub> receptor agonist. *J. Pharmacol. Exp. Ther.* **304(3)**, 1209-1216 (2003).

### 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, 11/04/2022

## 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