

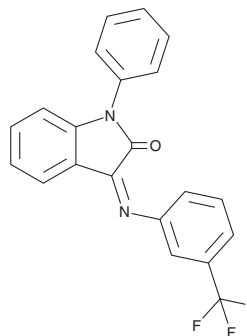
# PRODUCT INFORMATION



**HT-2157**

Item No. 33848

**CAS Registry No.:** 303149-14-6  
**Formal Name:** 1,3-dihydro-1-phenyl-3-[[3-(trifluoromethyl)phenyl]imino]-2H-indol-2-one  
**Synonym:** SNAP 37889  
**MF:** C<sub>21</sub>H<sub>13</sub>F<sub>3</sub>N<sub>2</sub>O  
**FW:** 366.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 250 nm  
**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

HT-2157 is supplied as a solid. A stock solution may be made by dissolving the HT-2157 in the solvent of choice, which should be purged with an inert gas. HT-2157 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of HT-2157 in ethanol is approximately 5 mg/ml and approximately 30 mg/ml in DMSO and DMF.

## Description

HT-2157 is an antagonist of the galanin-3 receptor (GAL<sub>3</sub>; K<sub>i</sub> = 17 nM).<sup>1</sup> It is selective for GAL<sub>3</sub> over GAL<sub>1</sub> and GAL<sub>2</sub> (K<sub>i</sub>s = >10 μM for both). HT-2157 (10 μM) induces apoptosis in various cell types, including HL-60 leukemia and BV-2 murine microglial cells, as well as isolated human peripheral blood mononuclear cells (PBMCs).<sup>2</sup> It decreases operant alcohol, sucrose, and saccharin self-administration in rats when administered at a dose of 30 mg/kg.<sup>3</sup> HT-2157 (3 and 10 mg/kg) increases drinking in the Vogel punished drinking task and reduces immobility time in the forced swim test in rats, indicating anxiolytic- and antidepressant-like activities, respectively.<sup>4</sup>

## References

1. Konkel, M.J., Lagu, B., Boteju, L.W., *et al.* 3-arylimino-2-indolones are potent and selective galanin GAL<sub>3</sub> receptor antagonists. *J. Med. Chem.* **49**(13), 3757-3758 (2006).
2. Koller, A., Rid, R., Beyreis, M., *et al.* *In vitro* toxicity of the galanin receptor 3 antagonist SNAP 37889. *Neuropeptides* **56**, 83-88 (2016).
3. Ash, B.L., Zanatta, S.D., Williams, S.J., *et al.* The galanin-3 receptor antagonist, SNAP 37889, reduces operant responding for ethanol in alcohol-preferring rats. *Regul. Pept.* **166**(1-3), 59-67 (2011).
4. Swanson, C.J., Blackburn, T.P., Zhang, X., *et al.* Anxiolytic- and antidepressant-like profiles of the galanin-3 receptor (Gal<sub>3</sub>) antagonists SNAP 37889 and SNAP 398299. *Proc. Nat. Acad. Sci. USA* **102**(48), 17489-17494 (2005).

### 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, 11/11/2022

## 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](http://WWW.CAYMANCHEM.COM)