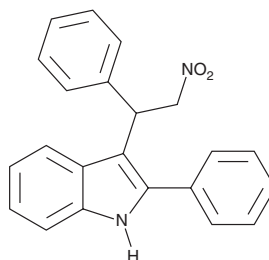


# PRODUCT INFORMATION

## GAT211

Item No. 24484

**CAS Registry No.:** 102704-40-5  
**Formal Name:** 3-(2-nitro-1-phenylethyl)-2-phenyl-1H-indole  
**MF:** C<sub>22</sub>H<sub>18</sub>N<sub>2</sub>O<sub>2</sub>  
**FW:** 342.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 300 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

GAT211 is supplied as a crystalline solid. A stock solution may be made by dissolving the GAT211 in the solvent of choice, which should be purged with an inert gas. GAT211 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of GAT211 in these solvents is approximately 5, 10, and 25 mg/ml, respectively.

GAT211 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GAT211 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. GAT211 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

GAT211 is an agonist and positive allosteric modulator (PAM) of cannabinoid receptor 1 (CB<sub>1</sub>) and a racemic mixture of GAT228 (Item No. 24485) and GAT229 (Item No. 24486), which have enantiomer-specific activities.<sup>1</sup> GAT211 increases β-arrestin recruitment and cAMP inhibition in HEK293A cells expressing GFP-tagged human CB<sub>1</sub> (hCB<sub>1</sub>-GFP) in a concentration-dependent manner. It also enhances the binding of the CB<sub>1</sub> full agonist CP 55,940 to CHO cells expressing hCB<sub>1</sub>, as well as the activity of 2-arachidonoyl glycerol (2-AG; Item No. 62160), arachidonoyl ethanolamide (AEA; Item No. 90050), and CP 55,940 in arrestin2 recruitment assays and increases ERK1/2 and PLCβ3 phosphorylation in HEK293 cells expressing hCB<sub>1</sub>. It inhibits excitatory postsynaptic currents (EPSCs) in a subset of CB<sub>1</sub>-expressing murine autaptic hippocampal neurons when used at a concentration of 1 μM.<sup>2</sup> GAT211 also decreases mechanical hypersensitivity in wild-type (EC<sub>50</sub> = 9.75 mg/kg), but not CB<sub>1</sub> knockout, mice in a model of inflammatory pain induced by complete Freund's adjuvant (CFA), and decreases mechanical and cold allodynia in a mouse model of paclitaxel-induced neuropathic pain when used at doses of 10 and 20 mg/kg per day.<sup>3</sup>

### References

1. Laprairie, R.B., Kulkarni, P.M., Deschamps, J.R., *et al.* Enantiospecific allosteric modulation of cannabinoid 1 receptor. *ACS Chem Biol.* **8**(6), 1188-1203 (2017).
2. Mitjavila, J., Yin, D., Kulkarni, P.M., *et al.* Enantiomer-specific positive allosteric modulation of CB<sub>1</sub> signaling in autaptic hippocampal neurons. *Pharmacol. Res.* **129**, 475-481 (2018).
3. Slivicki, R.A., Xu, Z., Kulkarni, P.M., *et al.* Positive allosteric modulation of CB1 suppresses pathological pain without producing tolerance or dependence. *Biol. Psychiatry* **84**(10), 722-733 (2018).

#### 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, 11/16/2022

#### 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