

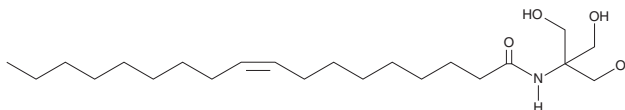
# PRODUCT INFORMATION



**JPC0323**

Item No. 39470

**CAS Registry No.:** 5972-45-2  
**Formal Name:** N-[2-hydroxy-1,1-bis(hydroxymethyl)ethyl]-9Z-octadecenamide  
**MF:** C<sub>22</sub>H<sub>43</sub>NO<sub>4</sub>  
**FW:** 385.6  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥2 years



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

## Laboratory Procedures

JPC0323 is supplied as a solid. A stock solution may be made by dissolving the JPC0323 in the solvent of choice, which should be purged with an inert gas. JPC0323 is soluble in chloroform and methanol.

## Description

JPC0323 is a positive allosteric modulator of the serotonin (5-HT) receptor subtypes 5-HT<sub>2A</sub> and 5-HT<sub>2C</sub>.<sup>1</sup> It is selective for these receptors over a panel of 21 receptors and transporters at 10 μM. JPC0323 (1 nM) enhances 5-HT-induced calcium release in CHO cells expressing recombinant human 5-HT<sub>2A</sub> or 5-HT<sub>2C</sub>. *In vivo*, JPC0323 (30 mg/kg) reduces spontaneous locomotor activity in rats.

## Reference

- Chen, J., Garcia, E.J., Merritt, C.R., *et al.* Discovery of novel oleamide analogues as brain-penetrant positive allosteric serotonin 5-HT<sub>2C</sub> receptor and dual 5-HT<sub>2C</sub>/5-HT<sub>2A</sub> receptor modulators. *J. Med. Chem.* **66**(14), 9992-10009 (2023).

### 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, 10/09/2023

## 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)