

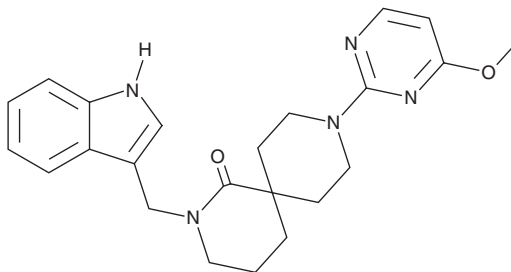
PRODUCT INFORMATION



IPSU

Item No. 30007

CAS Registry No.: 1373765-19-5
Formal Name: 2-(1H-indol-3-ylmethyl)-9-(4-methoxy-2-pyrimidinyl)-2,9-diazaspiro[5.5]undecan-1-one
MF: C₂₃H₂₇N₅O₂
FW: 405.5
Purity: ≥98%
UV/Vis.: λ_{max}: 220, 246 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

IPSU is supplied as a crystalline solid. A stock solution may be made by dissolving the IPSU in the solvent of choice, which should be purged with an inert gas. IPSU is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of IPSU in these solvents is approximately 1 mg/ml.

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

Description

IPSU is a potent orexin receptor 2 (OX2R) antagonist (K_i = 14.13 nM).¹ It is selective for OX2R over OX1R (K_i = 512.8 nM). IPSU (50 mg/kg) increases non-rapid eye movement (NREM) sleep in mice.

Reference

1. Betschart, C., Hintermann, S., Behnke, D., *et al.* Identification of a novel series of orexin receptor antagonists with a distinct effect on sleep architecture for the treatment of insomnia. *J. Med. Chem.* **56**(19), 7590-607 (2013).

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, 11/08/2022

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