

PRODUCT INFORMATION



PSB-KK1445

Item No. 42093

CAS Registry No.: 885896-54-8

Formal Name: 7-[(4-fluorophenyl)methyl]-3,7-dihydro-8-[[2-(1H-indol-3-yl)ethyl]amino]-1,3-dimethyl-1H-purine-2,6-dione

MF: $C_{24}H_{23}FN_6O_2$

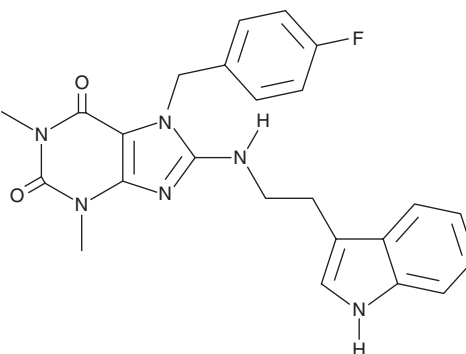
FW: 446.5

Purity: $\geq 95\%$

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

PSB-KK1445 is supplied as a solid. A stock solution may be made by dissolving the PSB-KK1445 in the solvent of choice, which should be purged with an inert gas. PSB-KK1445 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

PSB-KK1445 is an agonist of GPR18.¹ It is selective for GPR18 over GPR55 in β -arrestin recruitment assays using CHO-K1 cells expressing the human receptors ($EC_{50}s = 0.0454$ and $>10 \mu\text{M}$, respectively) and does not bind to cannabinoid 1 (CB_1) or CB_2 receptors ($K_i s = >10 \mu\text{M}$ for both).

Reference

1. Mahardhika, A.B., Załuski, M., Schoeder, C.T., *et al.* Potent, selective agonists for the cannabinoid-like orphan G protein-coupled receptor GPR18: A promising drug target for cancer and immunity. *J. Med. Chem.* **67**(12), 9896-9926 (2024).

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, 10/25/2024

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