

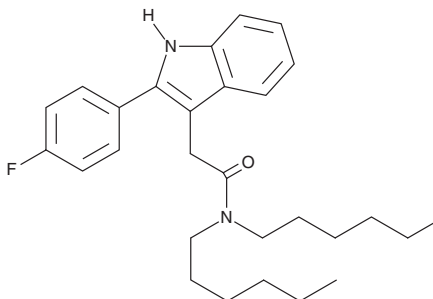
PRODUCT INFORMATION



FGIN-1-27

Item No. 18461

CAS Registry No.: 142720-24-9
Formal Name: 2-(4-fluorophenyl)-N,N-dihexyl-1H-indole-3-acetamide
MF: C₂₈H₃₇FN₂O
FW: 436.6
Purity: ≥98%
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 224, 301 nm
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

FGIN-1-27 is supplied as a crystalline solid. A stock solution may be made by dissolving the FGIN-1-27 in the solvent of choice. FGIN-1-27 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of FGIN-1-27 in ethanol and DMF is approximately 30 mg/ml and approximately 20 mg/ml in DMSO.

FGIN-1-27 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, FGIN-1-27 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. FGIN-1-27 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

FGIN-1-27 is an indoleacetamide that acts as a high affinity agonist of the 18 kDa translocator protein (TSPO; previously called the peripheral benzodiazepine receptor) with a K_i value of 5 nM.¹ It can enhance steroidogenesis of neuroactive steroids such as allopregnanolone (Item No. 16930).¹ Through this action, FGIN-1-27 indirectly potentiates GABA_A receptor signaling to produce anticonvulsant, anxiolytic, and sedative activity in both animal and clinical models.¹

Reference

1. James, M.L., Selleri, S., and Kassiou, M. Development of ligands for the peripheral benzodiazepine receptor. *Curr. Med. Chem.* **13**(17), 1991-2001 (2006).

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, 12/02/2022

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