

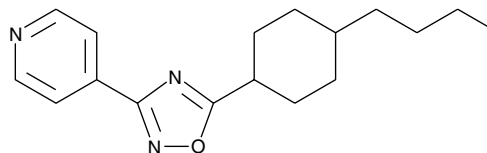
Product Information



PSN375963

Item No. 10008593

CAS Registry No.: 388575-52-8
Formal Name: 4-[5-(4-butylcyclohexyl)-1,2,4-oxadiazol-3-yl]-pyridine
MF: C₁₇H₂₃N₃O
FW: 285.4
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 224, 273 nm



Laboratory Procedures

For long term storage, we suggest that PSN375963 be stored as supplied at -20°C. It should be stable for at least two years.

PSN375963 is supplied as a crystalline solid. A stock solution may be made by dissolving the PSN375963 in an organic solvent purged with an inert gas. PSN375963 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of PSN375963 in ethanol and DMF is approximately 30 mg/ml and approximately 5 mg/ml in DMSO.

PSN375963 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, PSN375963 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. PSN375963 has a solubility of approximately 0.25 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

GPR119 (previously designated SNORF25) is an orphan G protein-coupled receptor expressed predominantly in the pancreas and gastrointestinal tract in humans and in the brain, pancreas, and gastrointestinal tract in rodents. It mediates a reduction in food intake and body weight gain in rats upon treatment with oleoyl ethanolamide (OEA), an endogenous, potent agonist for PPAR α .^{1,2} PSN375963 is a potent and selective agonist of GPR119 that shows similar potency to OEA at both recombinant mouse and human GPR119 receptors, exhibiting EC₅₀ values of 8.4 and 7.9 μ M, respectively (EC₅₀ values for OEA are 3.2 and 2.9 μ M, respectively).² These data suggest that PSN375963 may be useful as a therapeutic agent for the treatment of obesity.

References

1. Fu, J., Gaetani, S., Oveisi, F., *et al.* Oleylethanolamide regulates feeding and body weight through activation of the nuclear receptor PPAR- α . *Nature* **425**, 90-93 (2003).
2. Overton, H.A., Babbs, A.J., Doel, S.M., *et al.* Deorphanization of a G protein-coupled receptor for oleylethanolamide and its use in the discovery of small-molecule hypophagic agents. *Cell Metabolism* **3**, 167-175 (2006).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10008593

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY. NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog.**

Copyright Cayman Chemical Company, 02/14/2012

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com