

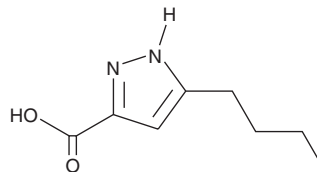
PRODUCT INFORMATION



LUF6283

Item No. 15423

CAS Registry No.: 92933-48-7
Formal Name: 5-butyl-1H-pyrazole-3-carboxylic acid
MF: C₈H₁₂N₂O₂
FW: 168.2
Purity: ≥98%
UV/Vis.: λ_{max}: 218 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

LUF6283 is supplied as a crystalline solid. A stock solution may be made by dissolving the LUF6283 in the solvent of choice, which should be purged with an inert gas. LUF6283 is soluble in organic solvents such as ethanol and DMSO. The solubility of LUF6283 in these solvents is approximately 16 mg/ml.

Description

LUF6283 is a partial agonist of GPR109A/HCA₂ G protein-coupled receptors (K_i = 0.55 μM in a radioligand binding assay).¹ It stimulates [³⁵S]-GTPγS binding and increases phosphorylation of ERK in HEK293T cells expressing GPR109A/HCA₂ receptors (EC₅₀s = 3.1 and 0.32 μM, respectively). *In vivo*, LUF6283 (400 mg/kg per day) reduces plasma levels of VLDL and hepatic expression of apolipoprotein B in mice.

Reference

1. Li, Z., Blad, C.C., van der Sluis, R.J., *et al.* Effects of pyrazole partial agonists on HCA₂-mediated flushing and VLDL-triglyceride levels in mice. *Br. J. Pharmacol.* **167**(4), 818-825 (2012).

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, 11/07/2022

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