

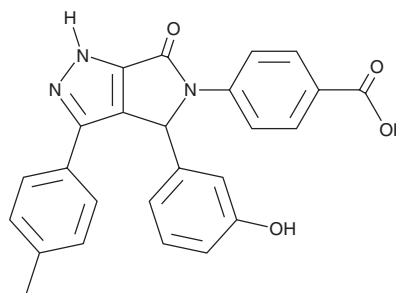
PRODUCT INFORMATION



CID-16020046

Item No. 15247

CAS Registry No.: 834903-43-4
Formal Name: 4-[4,6-dihydro-4-(3-hydroxyphenyl)-3-(4-methylphenyl)-6-oxopyrrolo[3,4-c]pyrazol-5(1H)-yl]-benzoic acid
MF: C₂₅H₁₉N₃O₄
FW: 425.4
Purity: ≥98%
UV/Vis.: λ_{max}: 280 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

CID-16020046 is supplied as a crystalline solid. A stock solution may be made by dissolving the CID-16020046 in the solvent of choice, which should be purged with an inert gas. CID-16020046 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CID-16020046 in these solvents is approximately 25 mg/ml.

Description

CID-16020046 is a GPR55 inverse agonist that antagonizes GPR55 constitutive activity with an IC₅₀ value of 15 μM.¹ It inhibits GPR55-mediated ERK1/2 phosphorylation, LPI-induced Ca²⁺ signaling (IC₅₀ = 0.21 μM in HEK-GPR55 cells), and GPR55-mediated transcription factor activation.¹ It does not affect ERK1/2 phosphorylation or transcription factor activation in CB receptor expressing cells and demonstrates weak activity against a broad spectrum of other GPCRs, ion channels, kinases, and nuclear receptors.¹ This compound has been shown to block GPR55-mediated endothelial wound healing and reverse LPI-inhibited platelet aggregation.¹

Reference

1. Kargl, J., Brown, A.J., Andersen, L., *et al.* A selective antagonist reveals a potential role of G protein-coupled receptor 55 in platelet and endothelial cell function. *J. Pharmacol. Exp. Ther.* **346**(1), 54-66 (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, 04/12/2023

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