

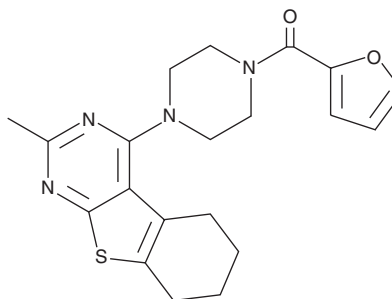
PRODUCT INFORMATION



ML-192

Item No. 15183

CAS Registry No.: 460331-61-7
Formal Name: 2-furanyl[4-(5,6,7,8-tetrahydro-2-methyl[1]benzothieno[2,3-d]pyrimidin-4-yl)-1-piperazinyl]-methanone
Synonym: CID-1434953
MF: C₂₀H₂₂N₄O₂S
FW: 382.5
Purity: ≥98%
UV/Vis.: λ_{max}: 235, 292 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

ML-192 is supplied as a crystalline solid. A stock solution may be made by dissolving the ML-192 in the solvent of choice, which should be purged with an inert gas. ML-192 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of ML-192 in these solvents is approximately 15, 5, and 10 mg/ml, respectively.

ML-192 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ML-192 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. ML-192 has a solubility of approximately 0.3 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.

Description

ML-192 is an antagonist of GPR55 (IC₅₀ = 702 nM).¹ It is selective for GPR55 over the cannabinoid (CB) receptors CB₁ and CB₂, as well as GPR35 (IC₅₀s = >32 μM).

Reference

1. Heynen-Genel, S., Dahl, R., Shi, S., *et al.* Selective GPR55 antagonists: Screening for selective ligands for GPR55. *Probe Reports from the NIH Molecular Libraries Program* (2011).

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, 10/12/2022

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