

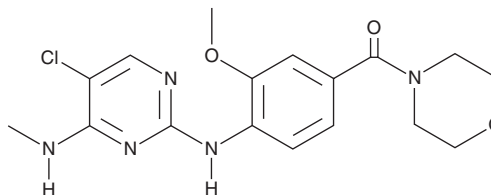
PRODUCT INFORMATION



HG-10-102-01

Item No. 20049

CAS Registry No.: 1351758-81-0
Formal Name: [4-[[5-chloro-4-(methylamino)-2-pyrimidinyl]amino]-3-methoxyphenyl]-4-morpholinyl-methanone
MF: C₁₇H₂₀ClN₅O₃
FW: 377.8
Purity: ≥98%
UV/Vis.: λ_{max}: 225, 275, 309 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

HG-10-102-01 is supplied as a crystalline solid. A stock solution may be made by dissolving the HG-10-102-01 in the solvent of choice. HG-10-102-01 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of HG-10-102-01 in these solvents is approximately 0.3, 30, and 10 mg/ml, respectively.

HG-10-102-01 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, HG-10-102-01 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. HG-10-102-01 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

HG-10-102-01 is an inhibitor of leucine-rich repeat kinase 2 (LRRK2; IC₅₀ = 20.3 nM).¹ It also inhibits the G2019S, A2016T, and [G2019S+A2016T] mutants of LRRK2 (IC₅₀s = 3.2, 153, and 95.9 nM, respectively).¹ HG-10-102-01 substantially inhibits Ser⁹¹⁰ and Ser⁹³⁵ phosphorylation of both wild-type and G2019S mutant LRRK2 in cells. When given intraperitoneally, HG-10-102-01 penetrates the blood-brain barrier and inhibits phosphorylation of LRRK2 in the brain as well as the kidney and spleen.¹

Reference

1. Choi, H.G., Zhang, J., Deng, X., *et al.* Brain Penetrant LRRK2 Inhibitor. *ACS Med. Chem. Lett.* **3(8)**, 658-662 (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, 12/20/2022

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