

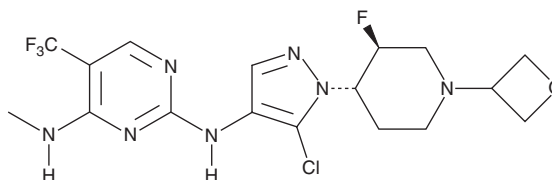
PRODUCT INFORMATION



GNE-9605

Item No. 23446

CAS Registry No.: 1536200-31-3
Formal Name: *rel*-N²-[5-chloro-1-[(3R,4R)-3-fluoro-1-(3-oxetanyl)-4-piperidinyl]-1H-pyrazol-4-yl]-N⁴-methyl-5-(trifluoromethyl)-2,4-pyrimidinediamine
MF: C₁₇H₂₀ClF₄N₇O
FW: 449.8
Purity: ≥98%
UV/Vis.: λ_{max}: 230 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

GNE-9605 is supplied as a crystalline solid. A stock solution may be made by dissolving the GNE-9605 in the solvent of choice. GNE-9605 is soluble in organic solvents such as methanol and DMSO, which should be purged with an inert gas. The solubility of GNE-9605 in these solvents is approximately 1 and 20 mg/ml, respectively.

Description

GNE-9605 is an orally bioavailable and potent inhibitor of leucine-rich repeat kinase 2 (LRRK2; IC₅₀ = 18.7 nM).¹ It is selective for LRRK2 over 178 kinases, inhibiting only TAK1-TAB1 >50% at a concentration of 0.1 μM. GNE-9605 (10 and 50 mg/kg) inhibits LRRK2 Ser¹²⁹² autophosphorylation in BAC transgenic mice expressing human LRRK2 protein with the G2019S mutation found in families with autosomal Parkinson's disease.

Reference

1. Estrada, A.A., Chan, B.K., Baker-Glenn, C., *et al.* Discovery of highly potent, selective, and brain-penetrant aminopyrazole leucine-rich repeat kinase 2 (LRRK2) small molecule inhibitors. *J. Med. Chem.* **57**(3), 921-936 (2014).

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, 10/17/2022

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