

PRODUCT INFORMATION

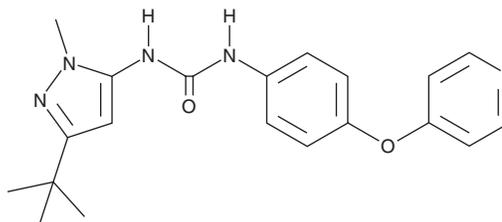


MAPK13 Inhibitor Compound 61

Item No. 42121

CAS Registry No.: 229002-10-2
Formal Name: N-[3-(1,1-dimethylethyl)-1-methyl-1H-pyrazol-5-yl]-N'-[4-(4-pyridinyloxy)phenyl]-urea
Synonyms: MAPK13-IN-1,
Mitogen-activated Protein Kinase 13 Inhibitor 1,
Mitogen-activated Protein Kinase 13 Inhibitor Compound 61

MF: C₂₀H₂₃N₅O₂
FW: 365.4
Purity: ≥98%
Supplied as: A 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

MAPK13 inhibitor compound 61 is supplied as a solid. A stock solution may be made by dissolving the MAPK13 inhibitor compound 61 in the solvent of choice, which should be purged with an inert gas. MAPK13 inhibitor compound 61 is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

MAPK13 inhibitor compound 61 is an inhibitor of p38 δ MAPK (IC₅₀ = 620 nM).¹

Reference

1. Yurtsever, Z., Patel, D.A., Kober, D.L., *et al.* First comprehensive structural and biophysical analysis of MAPK13 inhibitors targeting DFG-in and DFG-out binding modes. *Biochim. Biophys. Acta* **1860(11 Pt. A)**, 2335-2344 (2016).

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, 10/28/2024

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