

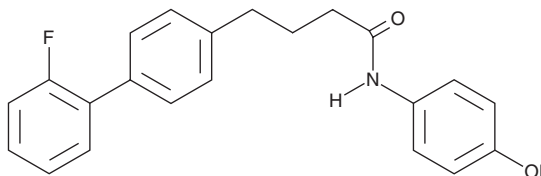
PRODUCT INFORMATION



MK2a Inhibitor

Item No. 17325

CAS Registry No.: 41179-33-3
Formal Name: 2'-fluoro-N-(4-hydroxyphenyl)-[1,1'-biphenyl]-4-butanamide
Synonym: CMPD 1
MF: C₂₂H₂₀FNO₂
FW: 349.4
Purity: ≥98%
UV/Vis.: λ_{max}: 249 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

MK2a inhibitor is supplied as a crystalline solid. A stock solution may be made by dissolving the MK2a inhibitor in the solvent of choice, which should be purged with an inert gas. MK2a inhibitor is soluble in organic solvents such as ethanol and DMSO. The solubility of MK2a inhibitor in these solvents is approximately 100 mM.

Description

MK2a inhibitor is a substrate-selective inhibitor of p38α MAPK that selectively inhibits p38α-dependent phosphorylation of MAPKAPK-2 (MK2) over activating transcription factor 2 (ATF2; K_i = 330 and >20,000 nM, respectively, in cell-free assays).¹ It inhibits proliferation of U87, A172, and U251 glioblastoma cells (EC₅₀s = 0.6-1 μM) *via* inhibition of tubulin polymerization and induction of apoptosis.²

References

1. Davidson, W., Frego, L., Peet, G.W., *et al.* Discovery and characterization of a substrate selective p38α inhibitor. *Biochemistry* **43**(37), 11658-11671 (2004).
2. Brennan, P.E. Deciphering the true antiproliferative target of an MK2 activation inhibitor in glioblastoma. *Cell Death Dis.* **e2069** (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, 02/17/2023

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