

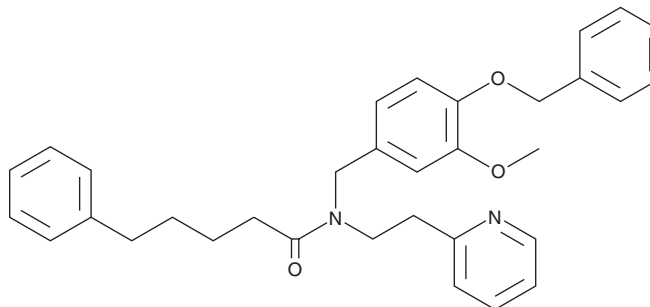
PRODUCT INFORMATION



MCP110

Item No. 37725

CAS Registry No.: 521310-51-0
Formal Name: N-[[3-methoxy-4-(phenylmethoxy)phenyl]methyl]-N-[2-(2-pyridinyl)ethyl]-benzenepentanamide
MF: C₃₃H₃₆N₂O₃
FW: 508.7
Purity: ≥95%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

MCP110 is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as acetonitrile and methanol purged with an inert gas can be used.

Description

MCP110 is an inhibitor of the protein-protein interaction between Ras and Raf-1.¹ It inhibits the interaction between H-Ras and Raf-1 in a yeast two-hybrid assay when used at a concentration of 30 μM. MCP110 (20 μM) inhibits Ras-induced activation of Raf-1 and MEK1 activity in HT-1080 fibrosarcoma cells. *In vivo*, MCP110 (100 and 600 mg/kg) reduces tumor volume in an SW620 colorectal cancer mouse xenograft model when administered alone or in combination with paclitaxel (Item No. 10461).²

References

1. Kato-Stankiewicz, J., Hakimi, I., Zhi, G., *et al.* Inhibitors of Ras/Raf-1 interaction identified by two-hybrid screening revert Ras-dependent transformation phenotypes in human cancer cells. *Proc. Natl. Acad. Sci. USA* **99(22)**, 14398-14403 (2002).
2. Skobeleva, N., Menon, S., Weber, L., *et al.* In vitro and in vivo synergy of MCP compounds with mitogen-activated protein kinase pathway- and microtubule-targeting inhibitors. *Mol. Cancer Ther.* **6(3)**, 898-906 (2007).

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, 12/13/2023

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