

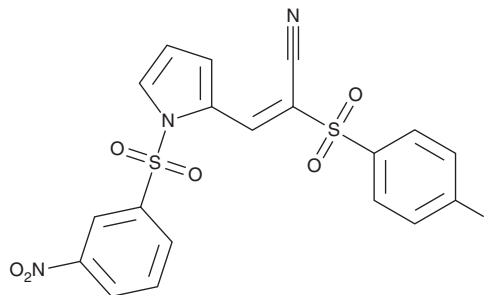
PRODUCT INFORMATION



AMZ30

Item No. 42052

CAS Registry No.: 1313613-09-0
Formal Name: (2E)-2-[(4-fluorophenyl)sulfonyl]-3-[1-[(3-nitrophenyl)sulfonyl]-1H-pyrrol-2-yl]-2-propenenitrile
Synonym: ML-136
MF: C₁₉H₁₂FN₃O₆S₂
FW: 461.4
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4years



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

Laboratory Procedures

AMZ30 is supplied as a crystalline solid. A stock solution may be made by dissolving the AMZ30 in the solvent of choice, which should be purged with an inert gas. AMZ30 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

AMZ30 is a covalent inhibitor of protein phosphatase methylesterase 1 (PME-1; IC₅₀ = 0.6 μM).¹ It reduces the levels of methylated protein phosphatase 2A (PP2A) in HEK293T cells stably overexpressing PME-1 when used at a concentration 20 μM. AMZ30 (210 μM) inhibits decreases in viability induced *tert*-butyl hydroperoxide (t-BHP) in U87MG and U-251MG glioblastoma cells.²

References

1. Bachovchin, D.A., Zuhl, A.M., Speers, A.E., *et al.* Discovery and optimization of sulfonyl acrylonitriles as selective, covalent inhibitors of protein phosphatase methylesterase-1. *J. Med. Chem.* **54**(14), 5229-5236 (2011).
2. Guffens, L., Derua, R., and Janssens, V. PME-1 sensitizes glioblastoma cells to oxidative stress-induced cell death by attenuating PP2A-B55α-mediated inactivation of MAPKAPK2-RIPK1 signaling. *Cell Death Discov.* **9**(1), 265 (2023).

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, 12/09/2024

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