

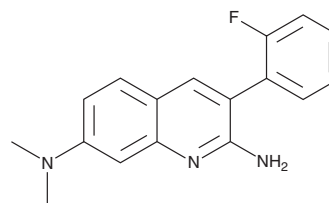
PRODUCT INFORMATION



Arylquin 1

Item No. 16961

CAS Registry No.: 1630743-73-5
Formal Name: 3-(2-fluorophenyl)-N⁷,N⁷-dimethyl-2,7-quinolinediamine
MF: C₁₇H₁₆FN₃
FW: 281.3
Purity: ≥98%
UV/Vis.: λ_{max}: 249, 267, 285 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

Arylquin 1 is supplied as a crystalline solid. A stock solution may be made by dissolving the arylquin 1 in the solvent of choice, which should be purged with an inert gas. Arylquin 1 is soluble in the organic solvent DMSO at a concentration of approximately 20 mg/ml.

Description

Arylquin 1 is a prostate apoptosis response-4 (Par-4) secretagogue.¹ It binds to vimentin to displace Par-4 for secretion. Arylquin 1 induces Par-4 secretion from mouse embryonic fibroblasts (MEFs) and prostate cancer cells *in vitro*, effects that can be blocked by the inhibitor of vesicular transport brefeldin A (Item No. 11861). Arylquin 1 induces paracrine apoptosis of H1299, HOP92, and H460 cancer cells when co-cultured with Par-4^{+/+}, but not Par-4^{-/-}, MEFs.

Reference

1. Burikhanov, R., Sviripa, V.M., Hebbar, N., *et al.* Arylquins target vimentin to trigger Par-4 secretion for tumor cell apoptosis. *Nat. Chem. Biol.* **10(11)**, 924-926 (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, 10/05/2022

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