

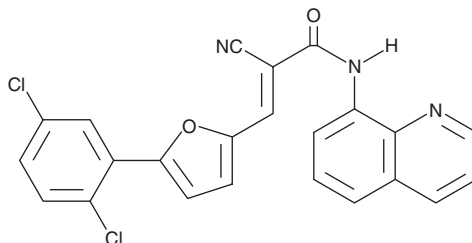
PRODUCT INFORMATION



AGK7

Item No. 16152

CAS Registry No.: 304896-21-7
Formal Name: 2-cyano-3-[5-(2,5-dichlorophenyl)-2-furanyl]-N-8-quinolinyl-2-propenamide
Synonym: SIRT2 Inhibitor (Inactive Control)
MF: C₂₃H₁₃Cl₂N₃O₂
FW: 434.3
Purity: ≥95%
UV/Vis.: λ_{max}: 219, 241, 391 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

AGK7 is supplied as a crystalline solid. A stock solution may be made by dissolving the AGK7 in the solvent of choice, which should be purged with an inert gas. AGK7 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of AGK7 in these solvents is approximately 0.5 and 0.2 mg/ml, respectively..

Description

AGK2 (Item No. 13145) is a cell-permeable, selective inhibitor of SIRT2 (IC₅₀ = 3.5 μM) that minimally affects either SIRT1 or SIRT3 at 10-fold higher levels.^{1,2} AGK7 is an inactive control to be used in experiments with AGK2. It differs from AGK2, structurally, only in the position of nitrogen in the quinoline group.¹ AGK7 has IC₅₀ values greater than 50 μM on SIRT1 and SIRT2 and greater than 5 μM on SIRT3.¹

References

1. Outeiro, T.F., Kontopoulos, E., Altmann, S.M., *et al.* Sirtuin 2 inhibitors rescue a-synuclein-mediated toxicity in models of Parkinson's Disease. *Science* **317(5837)**, 516-519 (2007).
2. Cole, P.A. Chemical probes for histone-modifying enzymes. *Nat. Chem. Biol.* **4(10)**, 590-597 (2008).

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, 11/22/2022

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