

PRODUCT INFORMATION



KYA1797K

Item No. 29221

CAS Registry No.: 1956356-56-1
Formal Name: (5Z)-5-[[5-(4-nitrophenyl)-2-furanylmethylene]-4-oxo-2-thioxo-3-thiazolidinepropanoic acid, monopotassium salt

MF: C₁₇H₁₁N₂O₆S₂ • K

FW: 442.5

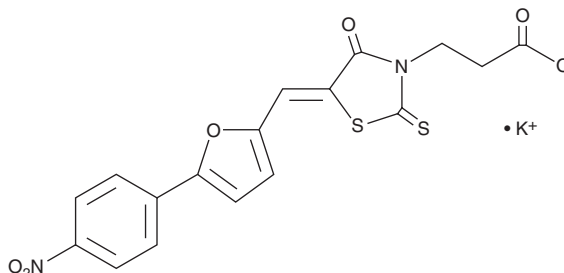
Purity: ≥95%

UV/Vis.: λ_{max}: 240, 327, 436 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

KYA1797K is supplied as a crystalline solid. A stock solution may be made by dissolving the KYA1797K in the solvent of choice, which should be purged with an inert gas. KYA1797K is soluble in the organic solvent DMSO at a concentration of approximately 0.5 mg/ml. KYA1797K is slightly soluble in ethanol and dimethyl formamide.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of KYA1797K can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of KYA1797K in PBS, pH 7.2, is approximately 0.3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

KYA1797K is an inhibitor of Wnt/β-catenin signaling.¹ It inhibits expression of a reporter induced by Wnt3a-conditioned medium in HEK293 cells (IC₅₀ = 0.75 μM). KYA1797K binds to the regulators of G-protein signaling (RGS) domain of axin and increases activation of GSK3β, as well as phosphorylation of β-catenin and Ras, in HEK293 cells when used at a concentration of 25 μM. It decreases β-catenin and Ras protein levels and inhibits colony formation in SW480 cells in a concentration-dependent manner. KYA1797K inhibits the growth of SW480, LoVo, DLD1, and HCT15 cells (GI₅₀s = 5, 4.8, 4.5, and 4.2 μM, respectively). It reduces tumor growth in a D-MT colorectal cancer mouse xenograft model when administered at a dose of 20 mg/kg.

Reference

1. Cha, P.-H., Cho, Y.-H., Lee, S.-K., *et al.* Small-molecule binding of the axin RGS domain promotes β-catenin and ras degradation. *Nat. Chem. Biol.* **12**(8), 593-600 (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, 10/21/2022

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