

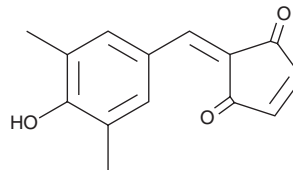
PRODUCT INFORMATION



TX-1918

Item No. 20840

CAS Registry No.: 503473-32-3
Formal Name: 2-[(4-hydroxy-3,5-dimethylphenyl)methylene]-4-cyclopentene-1,3-dione
MF: C₁₄H₁₂O₃
FW: 228.2
Purity: ≥98%
UV/Vis.: λ_{max}: 220, 248, 370 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

TX-1918 is supplied as a crystalline solid. A stock solution may be made by dissolving the TX-1918 in the solvent of choice, which should be purged with an inert gas. TX-1918 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). TX-1918 is slightly soluble in ethanol and DMF and soluble in DMSO at a concentration of approximately 0.33 mg/ml.

Description

TX-1918 is a cell-permeable tyrphostin derivative that inhibits eukaryotic elongation factor-2 kinase (eEF-2K; IC₅₀ = 440 nM).¹ It inhibits Src, PKA, PKC, and EGFR kinases at much higher concentrations (IC₅₀s = 4.4, 44, 44, and 440 μM, respectively).¹ TX-1918 disrupts the proliferation of HepG2 and HCT116 tumor cells (EC₅₀s = 2.07 and 230 μM, respectively) with ≥5,000-fold and ≥90-fold reduced mitochondrial- and hepato-toxicity, respectively when compared with the tyrphostin AG-17 (Item No. 10010248).¹

Reference

1. Hori, H., Nagasawa, H., Ishibashi, M., *et al.* TX-1123: An antitumor 2-hydroxyarylidene-4-cyclopentene-1,3-dione as a protein tyrosine kinase inhibitor having low mitochondrial toxicity. *Bioor. Med. Chem. Lett.* **10**(10), 3257-3265 (2002).

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, 09/30/2022

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