

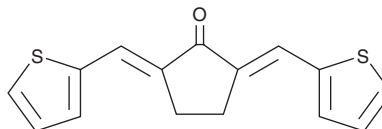
PRODUCT INFORMATION



CCT007093

Item No. 19992

CAS Registry No.: 176957-55-4
Formal Name: 2E,5E-bis(2-thienylmethylene)-cyclopentanone
MF: C₁₅H₁₂OS₂
FW: 272.4
Purity: ≥98%
UV/Vis.: λ_{max}: 255, 388 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

CCT007093 is supplied as a crystalline solid. A stock solution may be made by dissolving the CCT007093 in the solvent of choice, which should be purged with an inert gas. CCT007093 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of CCT007093 in these solvents is approximately 1 mg/ml.

CCT007093 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CCT007093 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CCT007093 has a solubility of approximately 0.33 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CCT007093 is an inhibitor of protein phosphatase 1D (PPM1D, also known as PP2Cδ and WIP1; IC₅₀ = 8.4 μM), which is encoded by a putative oncogene that is amplified in multiple cancer types.¹ It does not inhibit PPM1A. CCT007093 selectively inhibits the growth of cancer cells that overexpress PPM1D, including MCF-7, KPL-1, and MCF-3B cells, as well as certain ovarian clear cell carcinoma cells.^{1,2} CCT007093 has been used to elucidate the mechanism by which PPM1D drives growth in medulloblastoma cells.³

References

1. Rayter, S., Elliott, R., Travers, J., *et al.* A chemical inhibitor of PPM1D that selectively kills cells overexpressing PPM1D. *Oncogene* **27(8)**, 1036-1044 (2008).
2. Tan, D. S. P., Labmbros, M. B. K., Rayter, S., *et al.* PPM1D is a potential therapeutic target in ovarian clear cell carcinomas. *Clin. Cancer. Res* **15(7)**, 2269-2280 (2009).
3. Buss, M. C., Read, T.-A., Schniederjan, M. J., *et al.* HDM2 promotes WIP1-mediated medulloblastoma growth. *Neuro. Oncol.* **14(4)**, 440-458 (2012).

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, 10/25/2022

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