

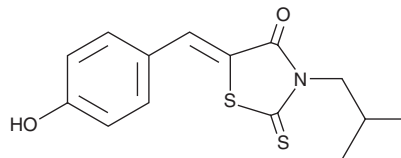
PRODUCT INFORMATION



PFM01

Item No. 36574

CAS Registry No.: 1558598-41-6
Formal Name: 5Z-[[4-hydroxyphenyl)methylene]-3-(2-methylpropyl)-2-thioxo-4-thiazolidinone
MF: C₁₄H₁₅NO₂S₂
FW: 293.4
Purity: ≥98%
Supplied as: A 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

PFM01 is supplied as a solid. A stock solution may be made by dissolving the PFM01 in the solvent of choice, which should be purged with an inert gas. PFM01 is soluble in the organic solvent DMSO at a concentration of approximately 29.3 mg/ml.

Description

PFM01 is an inhibitor of the endonuclease activity of the double-strand break repair protein Mre11.¹ It decreases the number of chromosomal breaks in 1BR3 fibroblasts induced by knockdown of BRCA2. PFM01 (50 μM) reduces homologous recombination-mediated DNA repair in DR-GFP U2OS cells and enhances DNA repair mediated by non-homologous end-joining (NHEJ) in H1299 dA3 cells.

Reference

1. Shibata, A., Moiani, D., Arvai, A.S., et al. DNA double-strand break repair pathway choice is directed by distinct MRE11 nuclease activities. *Mol. Cell* **53**(1), 7-18 (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, 11/23/2022

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