

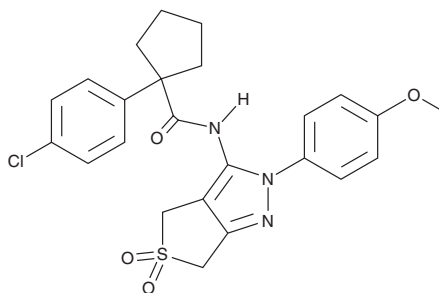
PRODUCT INFORMATION



TH301

Item No. 43814

CAS Registry No.: 450338-32-6
Formal Name: 1-(4-chlorophenyl)-N-[2,6-dihydro-2-(4-methoxyphenyl)-5,5-dioxido-4H-thieno[3,4-c]pyrazol-3-yl]-cyclopentanecarboxamide
MF: C₂₄H₂₄ClN₃O₄S
FW: 486.0
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

TH301 is supplied as a solid. A stock solution may be made by dissolving the TH301 in the solvent of choice, which should be purged with an inert gas. TH301 is slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO.

Description

TH301 is a stabilizer of cryptochrome circadian regulator 2 (CRY2).¹ It selectively stabilizes CRY2 over CRY1 in a degradation assay using HEK293 cells expressing CRY reporters at 0.1 to 10 μM. TH301 lengthens the circadian period in wild-type but not *Cry2*^{-/-} mouse fibroblasts. It enhances rosiglitazone-dependent upregulation of the brown adipose tissue-specific genes *Ucp1*, *Cidea*, and *Kcnk3* in primary mouse stromal vascular fractions.

Reference

1. Miller, S., Son, Y.L., Aikawa, Y., *et al.* Isoform-selective regulation of mammalian cryptochromes. *Nat. Chem. Biol.* **16**(6), 676-685 (2020).

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, 09/26/2025

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