

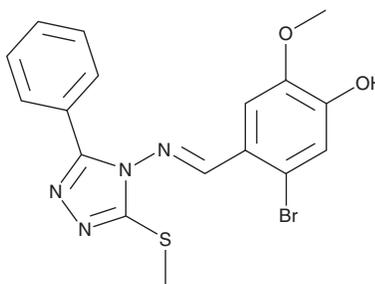
PRODUCT INFORMATION



GO289

Item No. 43959

CAS Registry No.: 694522-87-7
Formal Name: 5-bromo-2-methoxy-4-[[[3-(methylthio)-5-phenyl-4H-1,2,4-triazol-4-yl]imino]methyl]-phenol
MF: C₁₇H₁₅BrN₄O₂S
FW: 419.3
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

GO289 is supplied as a solid. A stock solution may be made by dissolving the GO289 in the solvent of choice, which should be purged with an inert gas. GO289 is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

GO289 is an inhibitor of casein kinase 2 (CK2; IC₅₀ = 7 nM).¹ It is selective for CK2 over a panel of 59 other kinases at 5 μM. GO289 (8 or 24 μM) increases the circadian period in a reporter assay using U2OS cells. It decreases levels of phosphorylated Akt and cell division cycle 37 (Cdc37) in C2C12 skeletal muscle myoblasts when used at concentrations of 1, 2, 3, or 4 μM.²

References

1. Oshima, T., Niwa, Y., Kuwata, K., *et al.* Cell-based screen identifies a new potent and highly selective CK2 inhibitor for modulation of circadian rhythms and cancer cell growth. *Sci. Adv.* **5**(1), eaau9060 (2019).
2. Borgo, C., Cesaro, L., Hirota, T., *et al.* Comparing the efficacy and selectivity of Ck2 inhibitors. A phosphoproteomics approach. *Eur. J. Med. Chem.* **214**, 113217 (2021).

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, 08/18/2025

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