

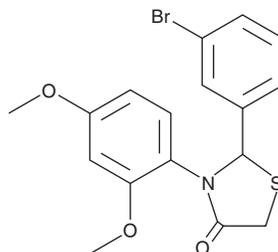
PRODUCT INFORMATION



CK-869

Item No. 15251

CAS Registry No.: 388592-44-7
Formal Name: 2-(3-bromophenyl)-3-(2,4-dimethoxyphenyl)-4-thiazolidinone
Synonym: CK-0157869
MF: C₁₇H₁₆BrNO₃S
FW: 394.3
Purity: ≥98%
UV/Vis.: λ_{max}: 278 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

CK-869 is supplied as a crystalline solid. A stock solution may be made by dissolving the CK-869 in the solvent of choice, which should be purged with an inert gas. CK-869 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CK-869 in these solvents is approximately 2.5 and 1.5 mg/ml, respectively.

CK-869 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CK-869 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CK-869 has a solubility of approximately 0.05 mg/ml in a 1:20 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CK-869 is an inhibitor of actin polymerization via inhibition of the actin-related protein 2/3 (Arp2/3) complex, leading to reduced cell motility rate and rearrangement of F-actin in M-1 cells.¹ It inhibits comet tail formation by *L. monocytogenes* with an IC₅₀ value of 7 μM.² It also impairs cell adhesion and cell spreading in MCF10A human breast epithelial cells.³

References

1. Ilatovskaya, D.V., Chubinskiy-Nadexhdin, V., Pavlov, T.S., *et al.* Arp2/3 inhibition induces amoeboid-like protrusions in MCF10A epithelial cells by reduced cytoskeletal-membrane coupling and focal adhesion assembly. *Cell Tissue Res.* **354**(3), 783-792 (2013).
2. Nolen, B.J., Tomasevic, N., Russell, A., *et al.* Characterization of two classes of small molecule inhibitors of Arp2/3 complex. *Nature* **460**(7258), 1031-1034 (2009).
3. Beckham, Y., Vasquez, R.J., Stricker, J., *et al.* Arp2/3 inhibition induces amoeboid-like protrusions in MCF10A epithelial cells by reduced cytoskeletal-membrane coupling and focal adhesion assembly. *PLoS One* **9**(6), e100943 (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, 11/16/2022

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