

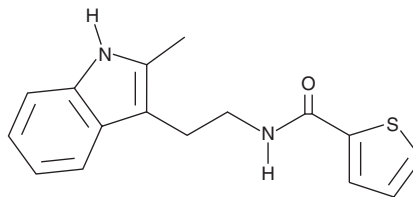
PRODUCT INFORMATION



CK-636

Item No. 16243

CAS Registry No.: 442632-72-6
Formal Name: N-[2-(2-methyl-1H-indol-3-yl)ethyl]-2-thiophenecarboxamide
MF: C₁₆H₁₆N₂OS
FW: 284.4
Purity: ≥98%
UV/Vis.: λ_{max}: 226, 273 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-636 is supplied as a crystalline solid. A stock solution may be made by dissolving the CK-636 in the solvent of choice, which should be purged with an inert gas. CK-636 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CK-636 in ethanol is approximately 20 mg/ml and approximately 30 mg/ml in DMSO and DMF.

CK-636 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CK-636 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CK-636 has a solubility of approximately 0.5 mg/ml in a 1:1 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-636 is an inhibitor of Arp2/3 complex action that binds the complex and inhibits actin polymerization *in vitro*, using human, bovine and fission yeast proteins (IC₅₀s = 4, 32, and 24 μM, respectively).¹ It is cell permeable, as it prevents actin polymerization and the formation of actin filament comet tails by *Listeria* in infected SKOV3 cells (IC₅₀ = 22 μM).¹ CK-636 is used to explore the role of Arp2/3 complex in cellular functions, including epidermal barrier formation and T cell migration.^{2,3}

References

1. 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).
2. Zhou, K., Muroyama, A., Underwood, J., *et al.* Actin-related protein2/3 complex regulates tight junctions and terminal differentiation to promote epidermal barrier formation. *Proc. Natl. Acad. Sci. USA* **110**(40), E3820-E3829 (2013).
3. Kwon, K.W., Park, H., and Doh, J. Migration of T cells on surfaces containing complex nanotopography. *PLoS One* **8**(9), e73960 (2013).

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, 11/28/2022

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