

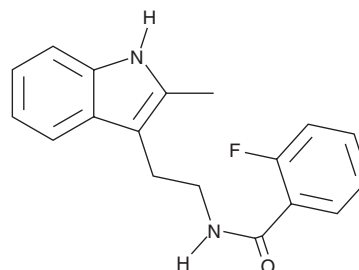
PRODUCT INFORMATION



CK-666

Item No. 29038

CAS Registry No.: 442633-00-3
Formal Name: 2-fluoro-N-[2-(2-methyl-1H-indol-3-yl)ethyl]-benzamide
MF: C₁₈H₁₇FN₂O
FW: 296.3
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-666 is supplied as a crystalline solid. A stock solution may be made by dissolving the CK-666 in the solvent of choice, which should be purged with an inert gas. CK-666 is soluble in ethanol and DMSO.

Description

CK-666 is an inhibitor of the actin-related protein 2/3 (Arp2/3) complex (IC₅₀s = 4, 17, and 5 μM for the human, bovine, and *S. pombe* complexes, respectively, in an actin polymerization assay).¹ It binds at the Arp2/3 interface, stabilizing the inactive conformation of the complex.² CK-666 (40 μM) inhibits actin comet tail formation around intracellular *L. monocytogenes* in infected SKOV3 cells.¹ It induces actin remodeling, inhibits lamellipodia formation, and reduces the motility rate in M-1 epithelial cells when used at a concentration of 200 μM.³ CK-666 (100 μM) impairs long-term fear memory formation in rats when microinjected into the lateral amygdala.⁴

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. Hetrick, B., Han, M.S., Helgeson, L.A., *et al.* Small molecules CK-666 and CK-869 inhibit actin-related protein 2/3 complex by blocking an activating conformational change. *Chem. Biol.* **20(5)**, 701-712 (2013).
3. Ilatovskaya, D.V., Chubinskiy-Nadexhdin, V., Pavlov, T.S., *et al.* Arp2/3 complex inhibitors adversely affect actin cytoskeleton remodeling in the cultured murine kidney collecting duct M-1 cells. *Cell Tissue Res.* **354(3)**, 783-792 (2013).
4. Basu, S., Kustanovich, I., and Lamprecht, R. Arp2/3 and VASP are essential for fear memory formation in lateral amygdala. *eNeuro* **3(6)**, pii:ENEURO.0302-16.2016 (2016).

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, 10/03/2022

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