

CK-666

Cell permeable inhibitor human ($IC_{50}=4\mu M$), fission yeast ($IC_{50}=5\mu M$), *S. cerevisiae* ($IC_{50}=12\mu M$) and bovine ($IC_{50}=17\mu M$) Arp2/3 complex between Arp2 and Arp3. Inhibits *Listeria* motility ($IC_{50}=7\mu M$).

Citations: 5

[View Online »](#)

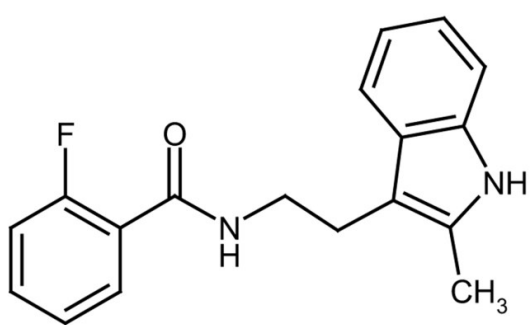
Ordering Information

[Order Online »](#)

ALX-270-506-M002	2mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Handling Protect from light.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 2-Fluoro-N-[2-(2-methyl-1H-indol-3-yl)ethyl]-benzamide

Appearance White to tan-yellow solid.

CAS 442633-00-3

Couple Target Arp

Couple Type Inhibitor

Formula $C_{18}H_{17}FN_2O$

MW 296.3

Purity ≥95% (HPLC)

Solubility Soluble in DMSO (a minimum concentration of 5 mg/ml is recommended).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)