

PD-98059

MEK inhibitor

PD-98059 is a potent and selective inhibitor of MAP kinase kinase (MEK). It selectively blocks the activation of MEK thereby inhibiting the phosphorylation and the activation of MAP kinase. PD-98059 is cell permeable. In PC12 pheochromocytoma cells, it completely blocked the increase in MAP kinase activity produced by NGF, $IC_{50}=2\mu M$. Enhances self-renewal of embryonic stem cells PD-98059 is an invaluable tool to help elucidate the role of the MAPK cascade in a variety of biological systems. Inhibits cell growth and proliferation in acute myelogenous leukemia cell lines by causing cell arrest.

Citations: 43

[View Online »](#)

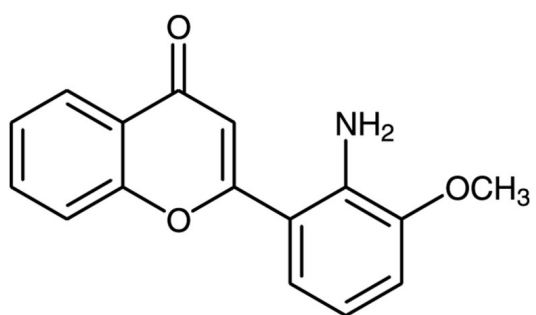
Ordering Information

[Order Online »](#)

BML-EI360-0005	5mg
----------------	-----

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

Appearance Pale yellow solid

CAS 167869-21-8

Couple Target MEK

Couple Type Inhibitor

Formula $C_{16}H_{13}NO_3$

MW 267.3

Purity ≥95% (HPLC)

Solubility Soluble in DMSO (6.5mg/ml); 25mg/ml in anhydrous DMSO), methanol, dichlormethane and 100% ethanol (0.6mg/ml).

Technical Info / Product Notes Replacement for **ADI-HPK-107**



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)