

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 »](#)

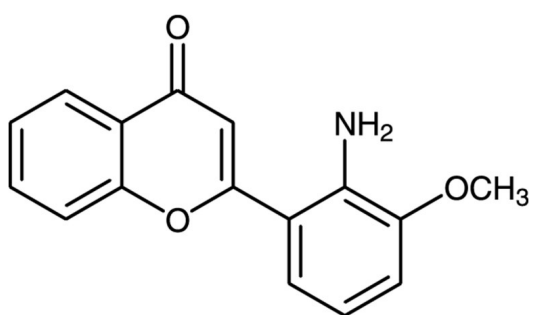
Ordering Information

[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**

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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