

U-0126

MEK inhibitor

U-0126 (Bis[amino[(2-aminophenyl)thio]methylene]butanedinitrile) is a novel, potent and selective MEK inhibitor (MEK1; IC_{50} =72nM, MEK2; IC_{50} =58nM). Inhibits AP-1 transcription activity. Blocks the production of inflammatory cytokines and matrix metalloproteinases.

Binding of extracellular growth factors to cell surface receptors often results in activation of the mitogen-activated protein kinase (MAPK). MAPK is regulated by MAPK kinase, also called MEK. Deprivation of growth factors during cell culture or intracellular MEK inhibition leads to inhibition of proliferation and apoptotic cell death. N Blank et al. (2002) have shown that U0126, a common used MEK inhibitor and compound that itself has no intrinsic fluorescence in vitro but develops an intensive fluorescence during cell culture which can be observed in all fluorescence channels with a predominance in the FL1 channel (525nm). Therefore either the use of PD98,059 or PD184352 as an alternative for U126 or nonfluorometric methods for detection of apoptosis should be considered.

Citations: 54

[View Online »](#)

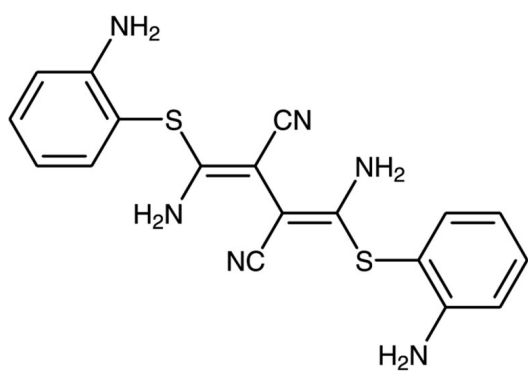
Ordering Information

[Order Online »](#)

BML-EI282-0001	1mg
BML-EI282-0005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at room temperature for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance	White solid
CAS	109511-58-2
Couple Target	MEK
Couple Type	Inhibitor
Formula	$C_{18}H_{16}N_6S_2$
MW	380.5
Purity	≥98% (TLC)
Purity Detail	In solution, U 0126 forms a mixture of isomers, but this does not affect biological activity.
Solubility	Soluble in DMSO (>10mg/ml) or 100% ethanol (2mg/ml).
Technical Info / Product Notes	Replacement for ADI-HPK-113



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