

SB239063

p38 inhibitor

A potent, cell permeable inhibitor of p38 MAP kinase (IC_{50} =44nM for recombinant purified human p38 α). Inhibits IL-1 and TNF- α production in LPS-stimulated human peripheral blood monocytes (IC_{50} =120nM and 350nM, respectively). Reported to protect against mild excitotoxic neuronal injury caused by NMDA and provides substantial protection against cell death induced by either oxygen glucose deprivation (OGD) or magnesium deprivation in cultured neurons. Also, reduces myocardial reperfusion injury via inhibition of endothelial adhesion molecule expression and blockade of polymorphonuclear (PMN) accumulation.

Citations: 9

[View Online »](#)

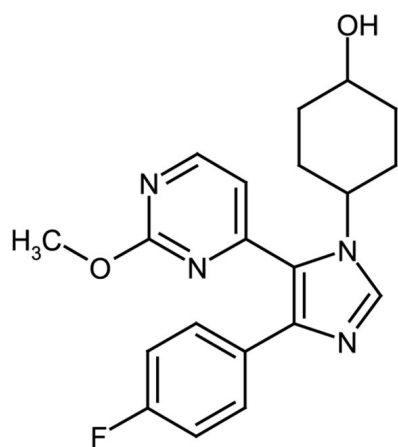
Ordering Information

[Order Online »](#)

ALX-270-351-M001	1mg
ALX-270-351-M005	5mg
ALX-270-351-C500	500 μ g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light. Packaged under inert gas.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	trans-1-(4-Hydroxycyclohexyl)-4-(fluorophenyl)-5-(2-methoxypyrimidin-4-yl) imidazole
Appearance	White to off-white solid.
CAS	193551-21-2
Couple Target	p38
Couple Type	Inhibitor
Formula	$C_{20}H_{21}FN_4O_2$
Identity	Identity determined by 1H -NMR.
MW	368.4
Purity	≥97% (HPLC)
Solubility	Soluble in DMSO, 100% ethanol or dichlormethane (may require slight warming).



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)