

Selumetinib

Potent and selective ATP non-competitive inhibitor of MEK1/2

Selumetinib (AZD6244) is a second-generation, orally available, potent and selective ATP non-competitive inhibitor of MEK 1/2. In direct assays, AZD6244 selectively inhibits purified active MEK1 and MEK2 with an IC₅₀ of 14 nM, and AZD6244 has reported IC₅₀ values of <100nM for MEK1 in ELISA assays and in bindings assays against recombinant domains of MEK2, Kd values were 53nM for AZD6244. AZD6244 demonstrated very effective anti-proliferation effects against several different cell lines with IC₅₀ values ranging from a low of 3nM to < 1 μ M depending upon the cell system used, and the compound inhibits basal and growth factor-stimulated phosphorylation of ERK1/ 2 with IC₅₀ concentrations < 40 nM. AZD6244 has demonstrated potent dose-dependent antitumor activity against a panel of mouse xenograft models of colorectal, pancreatic, liver, skin, and lung cancer, and inhibition of tumor growth was found to correlate with the reduction of phospho-ERK1/2 levels in tumors.

Citations: 1

[View Online »](#)

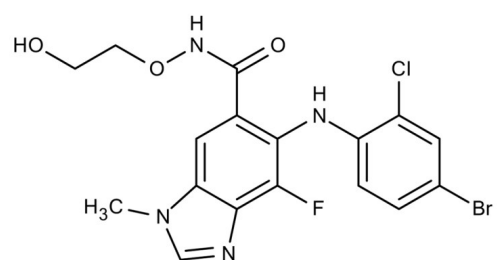
Ordering Information

[Order Online »](#)

ENZ-CHM184-0050	50mg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	AZD-6244, ARRY-142886
Appearance	Beige powder.
CAS	606143-52-6
Couple Target	MEK
Couple Type	Inhibitor
Formula	$C_{17}H_{15}BrClFN_4O_3$
Identity	Determined by EM-MS.
MW	457.7
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO (92mg/ml) or ethanol (<1mg/ml).



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)