

SMI-4a

Pim1 inhibitor

Selective inhibitor of Pim-1 and Pim-2 protein kinases. Inducer of G1 phase cell cycle arrest. Inducer of p27Kip1. Inducer of apoptosis through the mitochondrial pathway. Inhibitor of the mammalian target of rapamycin C1 (mTORC1) pathway. Downregulates c-myc. Inhibitor of PRAS40 phosphorylation and mTOR activity. Potential anti-cancer compound. Blocks prostate cancer growth.

Citations: 5

[View Online »](#)

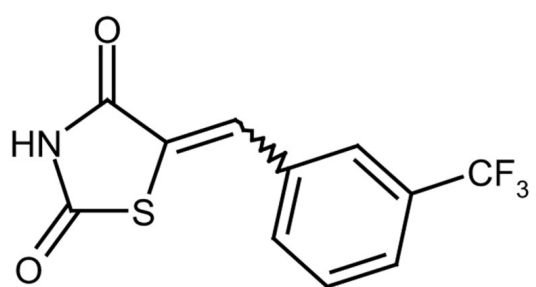
Ordering Information

[Order Online »](#)

ALX-270-515-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 6 months after receipt when stored at -20°C.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to off-white solid.
CAS	327033-36-3
Couple Target	Pim
Couple Type	Inhibitor
Formula	$C_{11}H_6F_3NO_2S$
MW	273.2
Purity	≥95% (HPLC, TLC, NMR/ELSD)
Solubility	Soluble in DMSO (at least 5mg/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-
fr@enzolifesciences.com](mailto:info-fr@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)