

AZD1208

pan-PIM kinase inhibitor

AZD1208 is a highly selective and potent pan-PIM kinase inhibitor. Displayed efficacy in models of acute myeloid leukemia, prostate cancer, and triple-negative breast cancer. Combination treatment of AZD1208 and imatinib eliminated chronic myelogenous leukemia stem cells in cell culture and a patient-derived mouse xenograft model. It improved immunotherapeutic antitumor T-cell response and disrupted the myeloid cell-mediated immunosuppressive tumor microenvironment increasing T-cell-mediated antitumor immunity.

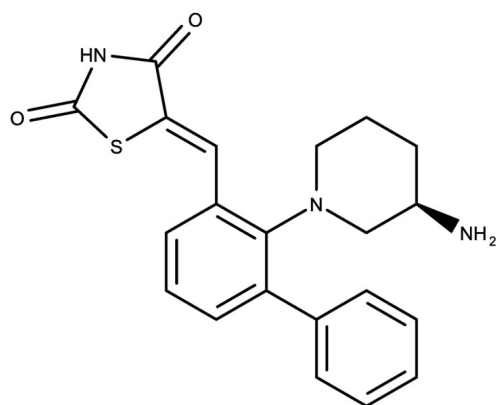
Ordering Information

[Order Online »](#)

ENZ-CHM457-0025	25mg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name (5Z)-5-[[2-[(3R)-3-Aminopiperidin-1-yl]-3-phenylphenyl]methylidene]-1,3-thiazolidine-2,4-dione

Appearance Yellow solid.

CAS 1204144-28-4

Couple Type Inhibitor

Formula $C_{21}H_{21}N_3O_2S$

Identity Determined by NMR.

MW 379.48

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (2 mg/ml with 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-
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)