

JNJ-28312141

Inhibitor of FMS

JNJ-28312141 is an orally active tyrosine kinase inhibitor for FMS with an IC_{50} of 0.69 nM and a cell assay IC_{50} of 2.6 nM.

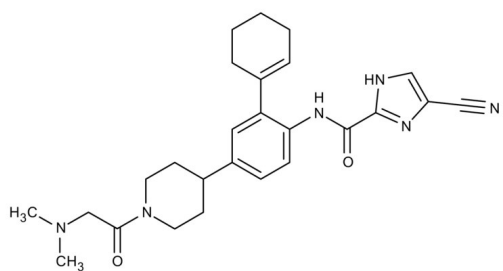
Ordering Information

[Order Online »](#)

ENZ-CHM190-0001	1mg
ENZ-CHM190-0005	5mg

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

Appearance	White powder.
CAS	885692-52-4
Couple Target	cFMS receptor tyrosine kinase
Couple Type	Inhibitor
Formula	$C_{26}H_{32}N_6O_2$
Identity	Determined by EM-MS, NMR
MW	460.6
Purity	≥99% (HPLC)
Solubility	Soluble in DMSO (12mg/ml) or ethanol (4mg/ml).

Last modified: May 29, 2024



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)