

Ruxolitinib

JAK inhibitor

Ruxolitinib is a Janus kinase inhibitor with IC_{50} values of 2.7 and 4.5, respectively, for JAK1 and JAK2.

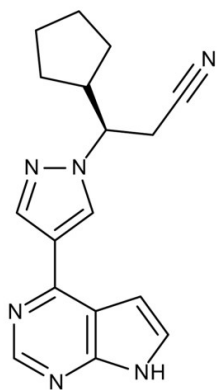
Ordering Information

[Order Online »](#)

ENZ-CHM156-0005	5mg
ENZ-CHM156-0010	10mg

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	INCB018424
Appearance	White powder.
CAS	941678-49-5
Couple Target	JAK
Couple Type	Inhibitor
Formula	$C_{17}H_{18}N_6$
Identity	Determined by EM-MS, NMR
MW	306.4
Purity	≥99% (HPLC)
Solubility	Soluble in DMSO (62mg/ml) or ethanol (2mg/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-
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)