

Baricitinib phosphate

Selective JAK1 and JAK2 inhibitor

Baricitinib is a selective JAK1 and JAK2 inhibitor. IC_{50} values are 5.9nM for JAK1, and 5.7nM for JAK2. In cell based assays Baricitinib/INCB-28050 demonstrated inhibition of intracellular signaling of multiple pro-inflammatory cytokines including IL-6 and IL-23 at concentrations of < 50 nM, and also showed good efficacy in a number of animal models at doses of 10mg/kg and below. Baricitinib is in Phase III development as a potential treatment for rheumatoid arthritis. It is in Phase II development as a potential treatment for psoriasis and diabetic nephropathy.

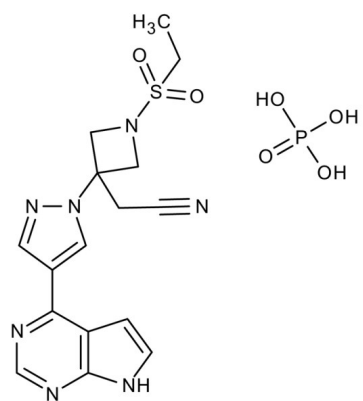
Ordering Information

[Order Online »](#)

ENZ-CHM198-0001	1mg
-----------------	-----

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	INCB028050, LY3009104, 2-[1-Ethylsulfonyl-3-[4-(7H-pyrrolo[2,3-d]pyrimidin-4-yl)pyrazol-1-yl]azetidin-3-yl]acetonitrile
Appearance	White powder.
CAS	1187595-84-1
Couple Target	JAK
Couple Type	Inhibitor
Formula	$C_{16}H_{17}N_7O_2S \cdot H_3PO_4$
Identity	Determined by EM-MS, NMR
MW	469.4
Purity	≥99% (HPLC)
Solubility	Soluble in DMSO (74mg/ml) or ethanol (1mg/ml).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com