

Product Information

Description:	INCB018424 (Ruxolitinib) is an orally bioavailable JAK inhibitor. (IC50 values for JAK1, JAK2 and JAK3 are 3, 5 and 332 nM, respectively). INCB018424 is useful for rheumatoid arthritis.
Chemical Name:	(3R)-3-cyclopentyl-3-[4-(7H-pyrrolo[2,3-d]pyrimidin-4-yl)pyrazol-1-yl]propanenitrile;phosphoric acid
Molecular Formula:	C ₁₇ H ₂₁ N ₆ O ₄ P
Supplied As:	Solid
Solubility:	≥20.2 mg/mL in DMSO; ≥6.92 mg/mL in EtOH with gentle warming and ultrasonic; ≥8.03 mg/mL in H ₂ O with gentle warming and ultrasonic
Stability:	At least 6 months at -20°C. Avoid freeze/thaw cycles.
Storage:	-20°C
MW:	404.36
Cas No.	1092939-17-7
Purity:	≥98% determined by HPLC

Chemical Structure

