

Ruxolitinib

JAK1 and JAK2 Inhibitor

Catalog code: tlr1-rux

<https://www.invivogen.com/ruxolitinib>

For research use only

Version 21J05-MM

PRODUCT INFORMATION

Contents

- 5 mg of Ruxolitinib

Storage and stability

- Ruxolitinib is shipped at room temperature. Upon receipt, store at -20°C.
- Upon resuspension in DMSO, prepare aliquots of ruxolitinib and store at -20°C. Avoid repeated freeze-thaw cycles. Resuspended product is stable for 6 months when properly stored.

Quality Control:

- Purity: ≥97% (UHPLC)
- The biological activity has been confirmed using cellular assays.
- The absence of bacterial contamination (e.g. lipoproteins and endotoxins) has been confirmed using HEK-Blue™ TLR2 and HEK-Blue™ TLR4 cells.

DESCRIPTION

Ruxolitinib (also known as INCB018424) is a potent, reversible, and selective Janus Kinase (JAK) 1 and JAK2 inhibitor¹. JAKs are constitutively bound to cytokine receptors, such as Type I interferons (IFNs) and interleukin-6 (IL-6). Upon binding of the ligand to the receptor, JAKs phosphorylate downstream targets such as STAT3/5, Akt, and ERK. This induces the production of cytokines and chemokines, including IFN-stimulated genes (ISGs). JAK-STAT signaling is crucial for the regulation and homeostasis of hematopoiesis and immunity². Ruxolitinib was developed as a potential therapeutic for a family of blood cancers termed myeloproliferative neoplasms (MPNs), which are characterized by the aberrant activation of the JAK-STAT pathway due to a mutation (V617F) in JAK2^{1,3}.

Ruxolitinib competes with ATP for binding to the catalytic site in the kinase domain, and thus inhibits not only the mutated JAK2 but wild-type JAK1-2 signaling pathways. Inhibition of the JAK-STAT signaling pathway by Ruxolitinib results in a dramatic decrease in levels of inflammatory cytokines, such as IL-6 and TNF-α. It is this attenuation of the inflammatory response that the clinical efficiency of Ruxolitinib is attributed². Ruxolitinib is approved for the treatment of the MPNs, myelofibrosis and polycythemia vera³. Pre-clinical data suggest that Ruxolitinib shows potential in the treatment of inflammatory conditions such as acute graft versus host disease⁴. Additionally, synergy has been reported between Ruxolitinib and the chemotherapy drug, dexamethasone, in the treatment of acute lymphoblastic leukemia in both *in vitro* and *in vivo* pre-clinical models⁵.

1. Quintas-Cardama, A. *et al.*, 2010. Preclinical characterization of the selective JAK1/2 inhibitor INCB018424: therapeutic implications for the treatment of myeloproliferative neoplasms. *Blood* 115, 3109-17. 2. Ajayi S. *et al.*, 2018. Ruxolitinib. *Recent Results Cancer Res* 212, 119-132. 3. Mascarenhas J. & Hoffman R. 2012. Ruxolitinib: the first FDA approved therapy for the treatment of myelofibrosis. *Clin Cancer Res* 18:3008-14. 4. Zeiser R. *et al.*, 2020. Ruxolitinib for Glucocorticoid-Refractory Acute Graft-versus-Host Disease. *N Engl J Med* 382:1800-10. 5. Verbeke D. *et al.*, 2019. Ruxolitinib Synergizes With Dexamethasone for the Treatment of T-cell Acute Lymphoblastic Leukemia. *Hemasphere* 3:e310.

CHEMICAL PROPERTIES

Structure:

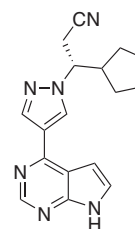
Synonym: INCB018424

Solubility: 10 mg/ml in DMSO

CAS number: 941678-49-5

Formula: C₁₇H₁₈N₆

Molecular weight: 306.37 g/mol



METHODS

Preparation of Ruxolitinib stock solution (20 mM)

1. Add 816 µl DMSO to 5 mg ruxolitinib and vortex until completely resuspended.
2. Prepare aliquots and store at -20°C. Prepare further dilutions using aqueous buffers. Do not store aqueous dilutions for more than 1 day.

Working concentration: 100 nM-10 µM for cell culture assays

Inhibition assay

Described below is a protocol to study the JAK-STAT pathway in the interferon (IFN) regulatory factor-inducible secreted embryonic alkaline phosphatase (SEAP) reporter B16 melanocytes, B16-Blue™ ISG cells.

1. Add 20 µl of ruxolitinib (final concentration 100 nM-10 µM) per well of a flat-bottom 96-well plate.
2. Add 160 µl (~75,000 cells) of a B16-Blue™ ISG cell suspension per well.
3. Incubate for 1 hour at 37°C in a 5% CO₂ incubator.
4. Add 20 µl of murine IFN-β (final concentration 100-1000 U/ml).
5. Incubate the plate at 37 °C in a 5% CO₂ incubator for 18-24 hours.
6. Monitor SEAP production using a SEAP detection assay, such as QUANTI-Blue™ Solution.

RELATED PRODUCTS

Product	Description	Cat. Code
3-Methyladenine	PI3K inhibitor	tlrl-3ma
Amlexanox	TBK1/IKKε inhibitor	inh-amx
B16-Blue™ ISG Cells	Reporter melanocytes	bb-ifnabg
QUANTI-Blue™ Solution	SEAP detection medium	rep-qbs
SB203580	p38 MAP kinase inhibitor	tlrl-sb20
U0126	MEK1 and MEK2 Inhibitor	tlrl-u0126

TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Asia: +852 3622-3480

E-mail: info@invivogen.com