

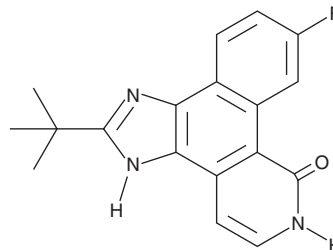
# PRODUCT INFORMATION



## JAK Inhibitor I

Item No. 15146

**CAS Registry No.:** 457081-03-7  
**Formal Name:** 2-(1,1-dimethylethyl)-9-fluoro-1,6-dihydro-7H-benz[h]imidazo[4,5-f]isoquinolin-7-one  
**Synonyms:** CMP 6, Janus-Associated Kinase Inhibitor I, Pyridone 6  
**MF:** C<sub>18</sub>H<sub>16</sub>FN<sub>3</sub>O  
**FW:** 309.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 219, 260, 270, 328, 343, 359 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

JAK inhibitor I is supplied as a crystalline solid. A stock solution may be made by dissolving the JAK inhibitor I in the solvent of choice, which should be purged with an inert gas. JAK inhibitor I is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of JAK inhibitor I in ethanol is approximately 1 mg/ml and approximately 15 mg/ml in DMSO and DMF.

JAK inhibitor I is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, JAK inhibitor I should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. JAK inhibitor I has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Janus-associated kinases (JAKs) are cytoplasmic tyrosine kinases that are required for activating the signaling of certain cytokines and growth factor receptors.<sup>1,2</sup> JAK1 mostly activates IL-6, whereas JAK1 and JAK3 trigger IL-2 and IL-4 and JAK1 and JAK2 stimulates IFN-γ. JAK Inhibitor I is a pyridine-containing tetracycle that interferes with JAK kinase activity by interacting within the ATP-binding cleft. It inhibits JAK1, 2, and 3 with IC<sub>50</sub> values of 15, 1, and 5 nM, respectively, while displaying significantly weaker affinities (IC<sub>50</sub>s = 130 nM - >10 μM) for other protein tyrosine kinases.<sup>3</sup> It was shown to block IL-2 and IL-4-dependent proliferation of mouse T-cell lymphoma cells with IC<sub>50</sub> values of 50-100 nM.<sup>3</sup> In a mouse model of atopic dermatitis, JAK Inhibitor I, supplied at 2 mg encapsulated in a PLGA nanoparticle, was shown to suppress IFN-γ/STAT1, IL-2/STAT5, and IL-4/STAT6 signaling pathways.<sup>4</sup>

### References

1. Leonard, W.J. and O'Shea, J.J. JAKS and STATS: Biological implications. *Annu. Rev. Immunol.* **16**, 293-322 (1998).
2. Yamaoka, K., Saharinen, P., Pesu, M., *et al.* The janus kinases (Jaks). *Genome Biol.* **5(12)**, 1-6 (2004).
3. Thompson, J.E., Cubbon, R.M., Cummings, R.T., *et al.* Photochemical preparation of a pyridone containing tetracycle: A Jak protein kinase inhibitor. *Bioorg. Med. Chem. Lett.* **12(8)**, 1219-1223 (2002).
4. Nakagawa, R., Yoshida, H., Asakawa, M., *et al.* Pyridone 6, a pan-JAK inhibitor, ameliorates allergic skin inflammation of NC/Nga mice *via* suppression of Th2 and enhancement of Th17. *J. Immunol.* **187(9)**, 4611-4620 (2011).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/19/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM