

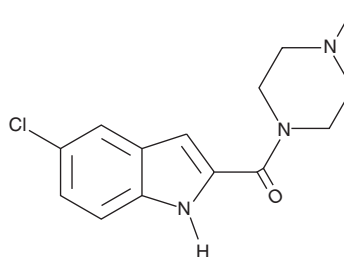
PRODUCT INFORMATION



JNJ-7777120

Item No. 10011925

CAS Registry No.: 459168-41-3
Formal Name: (5-chloro-1H-indol-2-yl)(4-methyl-1-piperazinyl)-methanone
MF: C₁₄H₁₆ClN₃O
FW: 277.8
Purity: ≥98%
UV/Vis.: λ_{max}: 219, 295 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

JNJ-7777120 is supplied as a crystalline solid. A stock solution may be made by dissolving the JNJ-7777120 in the solvent of choice. JNJ-7777120 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of JNJ-7777120 in these solvents is approximately 2 mg/ml in ethanol and approximately 14 mg/ml in DMSO and DMF.

JNJ-7777120 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, JNJ-7777120 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. JNJ-7777120 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

JNJ-7777120 is a potent and selective histamine H₄ receptor antagonist, with a K_i value of approximately 4 nM against the human, mouse, and rat H₄ receptors.¹ Its K_i values for the histamine H_{1,3} receptors exceed 1 μM, regardless of species, and it has no or negligible effects on a range of other receptors and transporters.¹ JNJ-7777120 inhibits mast cell chemotaxis induced by 10 μM histamine (IC₅₀ = 40 nM) and reduces neutrophil influx in mouse peritonitis models (10 mg/kg s.c.).¹ It also impairs eosinophil and lymphocyte influx into airways during allergic airway inflammation.^{2,3}

References

1. Thurmond, R.L., Desai, P.J., Dunford, P.J., *et al.* A potent and selective histamine H₄ receptor antagonist with anti-inflammatory properties. *J. Pharmacol. Exp. Ther.* **309**(1), 404-413 (2004).
2. Dunford, P.J., O'Donnell, N., Riley, J.P., *et al.* The histamine H₄ receptor mediates allergic airway inflammation by regulating the activation of CD4⁺ T cells. *J. Immunol.* **176**, 7062-7070 (2006).
3. Cowden, J.M., Riley, J.P., Ma, J.Y., *et al.* Histamine H₄ receptor antagonism diminishes existing airway inflammation and dysfunction via modulation of Th2 cytokines. *Respir. Res.* **11**(86), (2010).

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, 04/05/2024

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