

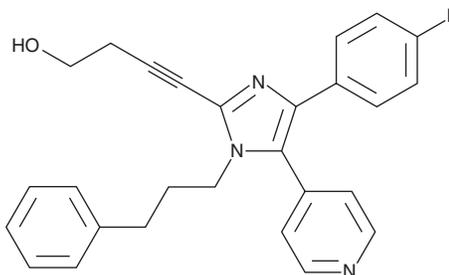
PRODUCT INFORMATION



RWJ 67657

Item No. 18718

CAS Registry No.: 215303-72-3
Formal Name: 4-[4-(4-fluorophenyl)-1-(3-phenylpropyl)-5-(4-pyridinyl)-1H-imidazol-2-yl]-3-butyn-1-ol
Synonym: JNJ-3026582
MF: C₂₇H₂₄FN₃O
FW: 425.5
Purity: ≥95%
UV/Vis.: λ_{max}: 241, 269 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

RWJ 67657 is supplied as a crystalline solid. A stock solution may be made by dissolving the RWJ 67657 in the solvent of choice. RWJ 67657 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of RWJ 67657 in these solvents is approximately 10, 5, and 2 mg/ml, respectively.

RWJ 67657 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, RWJ 67657 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. RWJ 67657 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

RWJ 67657 is an orally active inhibitor of the MAP kinases p38α and p38β (IC₅₀s = 1 and 11 μM, respectively, *in vitro*) that is inactive against p38γ and p38δ, as well as several other kinases.¹ It blocks the release of TNF-α and IL-1β from peripheral blood mononuclear cells stimulated with LPS (IC₅₀s = 3 and 11 nM, respectively) and inhibits TNF-α production in LPS-treated mice and rats.¹ RWJ 67657 potently blocks the proliferation of CD4⁺ peripheral blood T cells induced by CD28 stimulation alone (IC₅₀ = 0.5-4 nM).² It is commonly used to study the roles of p38α and p38β in cellular and whole animal systems.³⁻⁵

References

1. Wadsworth, S.A., Cavender, D.E., Beers, S.A., et al. *J. Pharmacol. Exp. Ther.* **291**(2), 680-687 (1999).
2. Schafer, P.H., Wadsworth, S.A., Wang, L., et al. *J. Immunol.* **162**(12), 7110-7119 (2016).
3. Kompa, A.R., See, F., Lewis, D.A., et al. *J. Pharmacol. Exp. Ther.* **325**(3), 741-750 (2008).
4. Antoon, J.W., Nitzchke, A.M., Martin, E.C., et al. *Int. J. Oncol.* **42**(4), 1139-1150 (2013).
5. Bai, Y.Y., Wang, L., Chang, D., et al. *Stroke* **46**(7), 1938-1946 (2015).

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, 11/08/2022

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