

PRODUCT INFORMATION



JNJ-0966

Item No. 31781

CAS Registry No.: 315705-75-0

Formal Name: N-[2-[(2-methoxyphenyl)amino]-4'-methyl[4,5'-bithiazol]-2'-yl]-acetamide

MF: C₁₆H₁₆N₄O₂S₂

FW: 360.5

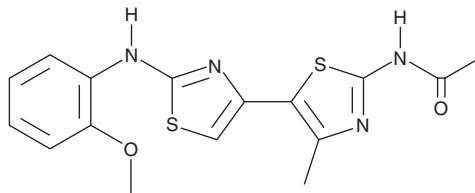
Purity: ≥98%

UV/Vis.: λ_{max}: 261, 311 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-0966 is supplied as a crystalline solid. A stock solution may be made by dissolving the JNJ-0966 in the solvent of choice, which should be purged with an inert gas. JNJ-0966 is soluble in methanol.

Description

JNJ-0966 is an inhibitor of matrix metalloproteinase-9 (MMP-9) zymogen activation (IC₅₀ = 440 nM).¹ It is selective for MMP-9 over MMP-1, MMP-2, and MMP-3 zymogen activation at 10 μM and does not inhibit the catalytic activity of MMP-1, -2, -3, -9, or -14. It inhibits invasion of HT-1080 cells in a Matrigel™ assay (IC₅₀ = 1 μM). JNJ-0966 (10 and 30 mg/kg, twice per day) reduces disease severity in a mouse model of experimental autoimmune encephalomyelitis (EAE).

Reference

1. Scannevin, R.H., Alexander, R., Haarlander, T.M., *et al.* Discovery of a highly selective chemical inhibitor of matrix metalloproteinase-9 (MMP-9) that allosterically inhibits zymogen activation. *J. Biol. Chem.* **292**(43), 17963-17974 (2017).

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, 11/20/2022

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