

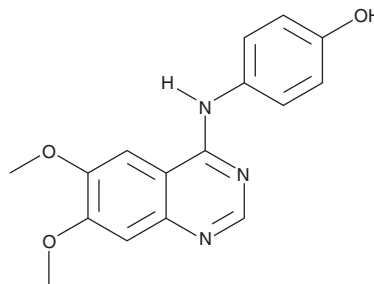
PRODUCT INFORMATION



WHI-P131

Item No. 10011246

CAS Registry No.: 202475-60-3
Formal Name: 4-[[6,7-dimethoxy-4-quinazoliny] amino]-phenol
Synonym: Janex 1
MF: C₁₆H₁₅N₃O₃
FW: 297.3
Purity: ≥98%
UV/Vis.: λ_{max}: 227, 247, 333 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

WHI-P131 is supplied as a crystalline solid. A stock solution may be made by dissolving the WHI-P131 in the solvent of choice. WHI-P131 is soluble in organic solvents such as methanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of WHI-P131 in these solvents is approximately 10, 1, and 1.5 mg/ml, respectively.

Description

Janus kinase 3 (JAK3) is a cytoplasmic protein tyrosine kinase that functions in the initiation of cytokine-triggered signal transduction by catalyzing the phosphorylation of signal transduction and activators of transcription (STATs). JAK3 is abundantly expressed in normal lymphoid cells as well as human lymphoblastic and myeloblastic leukemia cells. WHI-P131 selectively inhibits the tyrosine kinase activity of JAK3 with an IC₅₀ value of 78 μM without affecting the enzymatic activity of JAK1, JAK2, or other protein tyrosine kinases (IC₅₀ ≥350 μM).¹ WHI-P131 induces apoptosis in JAK3-expressing human leukemia cell lines NALM-6 and LC1;19, but not in melanoma or squamous carcinoma cell lines.¹

Reference

1. Sudbeck, E.A., Liu, X.-P., Narla, R.K., *et al.* Structure-based design of specific inhibitors of Janus kinase 3 as apoptosis-inducing antileukemic agents. *Clin. Cancer Res.* **5(6)**, 1569-1582 (1999).

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, 11/09/2022

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