

PRODUCT INFORMATION



CHIR98024

Item No. 34872

CAS Registry No.: 556813-39-9

Formal Name: N⁶-[2-[4-(2,4-dichlorophenyl)-5-(1H-imidazol-2-yl)-2-pyrimidinyl]amino]ethyl]-3-nitro-2,6-pyridinediamine

Synonym: CT 98024

MF: C₂₀H₁₇Cl₂N₉O₂

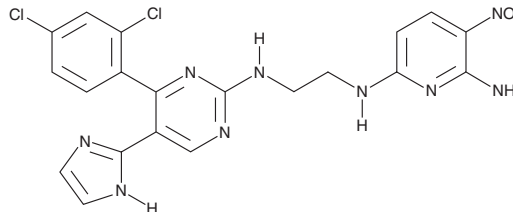
FW: 486.3

Purity: ≥95%

Supplied as: A 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

CHIR98024 is supplied as a solid. A stock solution may be made by dissolving the CHIR98024 in the solvent of choice, which should be purged with an inert gas. CHIR98024 is slightly soluble in ethanol and DMSO.

Description

CHIR98024 is an inhibitor of glycogen synthase kinase 3 (GSK3; IC₅₀ = 0.56 nM).¹

Reference

1. Wagman, A.S., Johnson, K.W., and Bussiere, D.E. Discovery and development of GSK3 inhibitors for the treatment of type 2 diabetes. *Curr. Pharm. Des.* **10(10)**, 1105-1137 (2004).

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, 12/08/2022

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