

PRODUCT INFORMATION

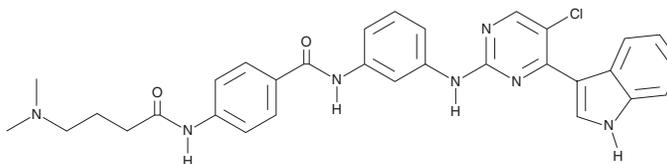


THZ1-R

Item No. 31054

CAS Registry No.: 1621523-07-6
Formal Name: N-[3-[[5-chloro-4-(1H-indol-3-yl)-2-pyrimidinyl]amino]phenyl]-4-[[4-(dimethylamino)-1-oxobutyl]amino]-benzamide

MF: C₃₁H₃₀ClN₇O₂
FW: 568.1
Purity: ≥98%
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

THZ1-R is supplied as a solid. A stock solution may be made by dissolving the THZ1-R in the solvent of choice, which should be purged with an inert gas. THZ1-R is soluble in the organic solvent DMSO at a concentration of approximately 10 mM.

Description

THZ1-R is a non-cysteine reactive derivative of the cyclin-dependent kinase 7 (Cdk7) inhibitor THZ1 (Item No. 9002215).¹ THZ1-R inhibits Cdk7 (IC₅₀s = 142-241 nM) less potently than THZ1 and lacks antiproliferative activity against T cell acute lymphoblastic leukemia (T-ALL) cell lines.

Reference

1. Kwiatkowski, N., Zhang, T., Rahl, P.B., *et al.* Targeting transcription regulation in cancer with a covalent CDK7 inhibitor. *Nature* **511**(7511), 616-620 (2014).

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, 12/09/2022

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