

PRODUCT INFORMATION



(R)-RP 6306

Item No. 41738

CAS Registry No.: 2719793-91-4
Formal Name: (1R)-2-amino-1-(3-hydroxy-2,6-dimethylphenyl)-5,6-dimethyl-1H-pyrrolo[2,3-b]pyridine-3-carboxamide

Synonym: (R)-Lunresertib

MF: C₁₈H₂₀N₄O₂

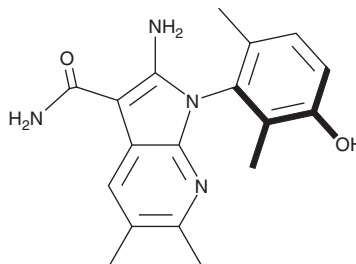
FW: 324.4

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

(R)-RP 6306 is supplied as a solid. A stock solution may be made by dissolving the (R)-RP 6306 in the solvent of choice, which should be purged with an inert gas. (R)-RP 6306 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

(R)-RP 6306 is a stereoisomer of RP 6306, an inhibitor of membrane-associated tyrosine- and threonine-specific Cdc2-inhibitory kinase (Myt1 kinase), also known as PKMYT1.¹ It inhibits Myt1 kinase with an IC₅₀ value of 1,360 nM.

Reference

1. Szychowski, J., Liu, B., Dietrich, E., *et al.* Compounds, pharmaceutical compositions, and methods of preparing compounds and their use. *Repare Therapeutics Inc.* US 2023/0122909 A1 (2023).

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, 09/16/2024

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