

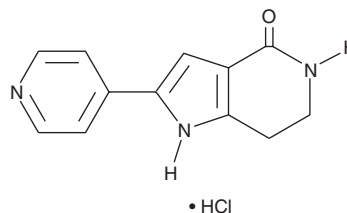
PRODUCT INFORMATION



PHA-767491 (hydrochloride)

Item No. 18218

CAS Registry No.: 942425-68-5
Formal Name: 1,5,6,7-tetrahydro-2-(4-pyridinyl)-4H-pyrrolo[3,2-c]pyridin-4-one, monohydrochloride
Synonym: CAY10572
MF: C₁₂H₁₁N₃O • HCl
FW: 249.7
Purity: ≥98%
UV/Vis.: λ_{max}: 216, 315 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

PHA-767491 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the PHA-767491 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. PHA-767491 (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of PHA-767491 (hydrochloride) in these solvents is approximately 5 and 0.3 mg/ml, respectively.

Description

PHA-767491 is a potent inhibitor of Cdc7 kinase with an IC₅₀ value of 10 nM in the presence of 1.5 μM ATP.¹ At 10 μM, PHA-767491 induces apoptotic cell death in multiple cancer cell types.¹ PHA-767491 also inhibits Cdk9, a kinase involved in the phosphorylation of RNA polymerase II and in transcriptional regulation of gene expression, with an IC₅₀ value of 34 nM.²

References

1. Montagnoli, A., Valsasina, B., Croci, V., *et al.* A Cdc7 kinase inhibitor restricts initiation of DNA replication and has antitumor activity. *Nat. Chem. Biol.* **4(6)**, 357-365 (2008).
2. Menichincheri, M., Bargiotti, A., Berthelsen, J., *et al.* First Cdc7 kinase inhibitors: Pyrrolopyridinones as potent and orally active antitumor agents. 2. lead discovery. *J. Med. Chem.* **52(2)**, 293-307 (2009).

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, 06/12/2023

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