

PRODUCT INFORMATION



ICEC0942 (hydrochloride)

Item No. 30997

CAS Registry No.: 1805789-54-1
Formal Name: (3R,4R)-4-[[[3-(1-methylethyl)-7-[(phenylmethyl)amino]pyrazolo[1,5-a]pyrimidin-5-yl]amino]methyl]-3-piperidinol, monohydrochloride

MF: C₂₂H₃₀N₆O • HCl

FW: 431.0

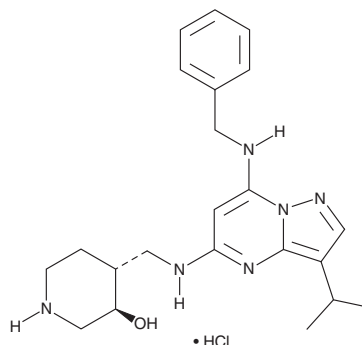
Purity: ≥95%

UV/Vis.: λ_{max}: 245 nm

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

ICEC0942 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the ICEC0942 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. ICEC0942 (hydrochloride) is soluble in the organic solvent DMSO.

Description

ICEC0942 is a cyclin-dependent kinase 7 (Cdk7) inhibitor (IC₅₀ = 0.04 μM).¹ It is selective for Cdk7 over Cdk1, -2, -4, -5, -6, and -9 (IC₅₀s = 1.8, 0.62, 49, 9.4, 34, and 1.2 μM, respectively). ICEC0942 induces cell cycle arrest at the G₂/M phase and apoptosis in HCT116 cells when used at concentrations of 1 and 10 μM. *In vivo*, ICEC0942 (100 mg/kg) reduces tumor growth in MCF-7 and HCT116 mouse xenograft models.

Reference

1. Patel, H., Periyasamy, M., Sava, G.P., *et al.* ICEC0942, an orally bioavailable selective inhibitor of CDK7 for cancer treatment. *Mol. Cancer Ther.* **17(6)**, 1156-1166 (2018).

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, 02/27/2023

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