

PRODUCT INFORMATION



SCR-1481B1

Item No. 33269

CAS Registry No.: 1174161-86-4
Formal Name: N-[4-[(2-amino-3-chloro-4-pyridinyl)oxy]-3-fluorophenyl]-5-(4-fluorophenyl)-4-oxo-1-[(phosphonoxy)methyl]-3(4H)-pyridinecarboxamide, compd. with 2-amino-2-(hydroxymethyl)-1,3-propanediol (1:2)

MF: C₂₄H₁₈ClF₂N₄O₇P • 2C₄H₁₁NO₃

FW: 821.1

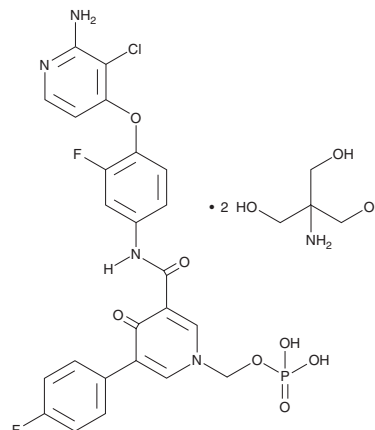
Purity: ≥95%

UV/Vis.: λ_{max}: 265 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

SCR-1481B1 is supplied as a crystalline solid. A stock solution may be made by dissolving the SCR-1481B1 in the solvent of choice, which should be purged with an inert gas. SCR-1481B1 is soluble in the organic solvent DMSO at a concentration of approximately 1 mg/ml.

SCR-1481B1 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SCR-1481B1 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. SCR-1481B1 has a solubility of approximately 0.11 mg/ml in a 1:8 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

SCR-1481B1 is a pyridone.¹

Reference

1. Furth, J.J. and Cohen, S.S. Inhibition of mammalian DNA polymerase by the 5'-triphosphate of 1-β-D-arabinofuranosylcytosine and the 5'-triphosphate of 9-β-D-arabinofuranoxyladenine. *Cancer Res.* **28(10)**, 2061-2067 (1968).

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, 11/14/2022

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