

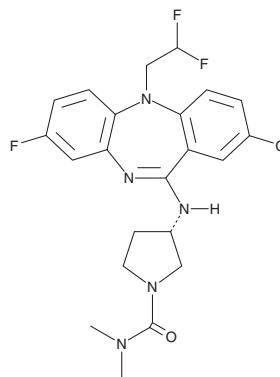
PRODUCT INFORMATION



NVS-PAK1-C

Item No. 19965

CAS Registry No.: 2250019-95-3
Formal Name: (3S)-(3-[[2-chloro-5-(2,2-difluoroethyl)-8-fluoro-5H-dibenzo[b,e][1,4]diazepin-11-yl]amino]-N,N-dimethyl-1-pyrrolidinecarboxamide
MF: C₂₂H₂₃ClF₃N₅O
FW: 465.9
Purity: ≥98%
UV/Vis.: λ_{max}: 210, 260, 298 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

NVS-PAK1-C is supplied as a crystalline solid. A stock solution may be made by dissolving the NVS-PAK1-C in the solvent of choice, which should be purged with an inert gas. NVS-PAK1-C is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of NVS-PAK1-C in ethanol is approximately 20 mg/ml and approximately 50 mg/ml in DMSO and DMF.

NVS-PAK1-C is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, NVS-PAK1-C should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. NVS-PAK1-C has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

NVS-PAK1-C is a negative control for the allosteric PAK1 inhibitor NVS-PAK1-1 (Item No. 19964). See the Structural Genomic Consortium (SGC) website for more information.

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, 10/25/2022

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