

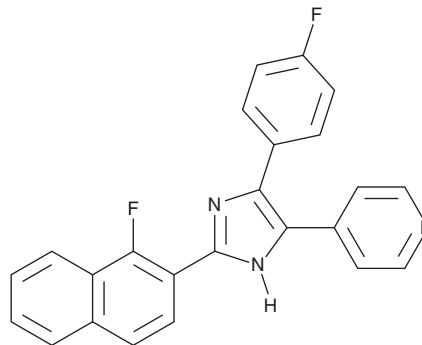
PRODUCT INFORMATION



CK1-IN-1

Item No. 37313

CAS Registry No.: 1784751-20-7
Formal Name: 4-[2-(1-fluoro-2-naphthalenyl)-4-(4-fluorophenyl)-1H-imidazol-5-yl]-pyridine
Synonyms: Casein Kinase 1 Inhibitor 1, CK1 Inhibitor 1
MF: C₂₄H₁₅F₂N₃
FW: 383.4
Purity: ≥98%
UV/Vis.: λ_{max}: 226, 279, 317 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

CK1-IN-1 is supplied as a solid. A stock solution may be made by dissolving the CK1-IN-1 in the solvent of choice, which should be purged with an inert gas. CK1-IN-1 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CK1-IN-1 in these solvents is approximately 1 and 3 mg/ml, respectively.

Description

CK1-IN-1 is an inhibitor of casein kinase 1δ (CK1δ) and CK1ε (IC₅₀s = 15 and 16 nM, respectively).¹ It is selective for CK1δ and CK1ε over TGF-β receptor type 1 (ALK5; IC₅₀ = 151.8 nM) and TGF-β receptor type 2 (TGFBR2) at 1 μM.

Reference

1. Zhong, Q., Laco, F., Liao, M.-C., *et al.* Influencing the fate of cardiac and neural stem cell differentiation using small molecule inhibitors of ALK5. *Stem Cells Transl. Med.* **7(10)**, 709-720 (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, 04/12/2023

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