

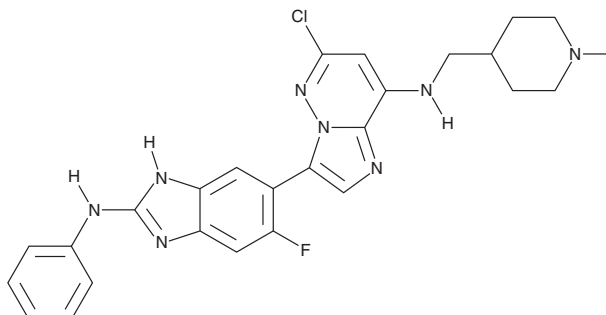
PRODUCT INFORMATION



IRE1 α Kinase-IN-1

Item No. 43728

CAS Registry No.: 2328097-41-0
Formal Name: 6-chloro-3-[6-fluoro-2-(phenylamino)-1H-benzimidazol-5-yl]-N-[(1-methyl-4-piperidinyl)methyl]-imidazo[1,2-b]pyridazin-8-amine
Synonyms: Inositol-requiring Enzyme 1 α Kinase-IN-1, Inositol-requiring Enzyme 1 α Kinase Inhibitor 1, IRE1 α Kinase Inhibitor 1
MF: C₂₆H₂₆ClFN₈
FW: 505.0
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

IRE1 α kinase-IN-1 is supplied as a solid. A stock solution may be made by dissolving the IRE1 α kinase-IN-1 in the solvent of choice, which should be purged with an inert gas. IRE1 α kinase-IN-1 is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

IRE1 α kinase-IN-1 is an allosteric inhibitor of inositol-requiring enzyme 1 α (IRE1 α ; IC₅₀s = 160 and 80 nM for IRE1 α kinase and ribonuclease activities, respectively).¹ It is selective for IRE1 α over a panel of 454 wild-type and mutant kinases at 1 μ M. IRE1 α kinase-IN-1 inhibits endoplasmic reticulum (ER) stress-induced IRE1 α oligomerization in HEK293 cells (IC₅₀ = 740 nM).

Reference

1. Colombano, G., Caldwell, J.J., Matthews, T.P., *et al.* Binding to an unusual inactive kinase conformation by highly selective inhibitors of inositol-requiring enzyme 1 α kinase-endoribonuclease. *J. Med. Chem.* **62**(5), 2447-2465 (2019).

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, 07/24/2025

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