

PRODUCT INFORMATION



FT709

Item No. 39971

CAS Registry No.: 2413991-74-7
Formal Name: (2S)-1-[5-[(2,3-dihydro-1,4-dioxino[2,3-b]pyridin-7-yl)sulfonyl]-3,4,5,6-tetrahydropyrrolo[3,4-c]pyrrol-2(1H)-yl]-2-hydroxy-2-(2-methyl-4-benzoxazolyl)-ethanone

MF: C₂₃H₂₂N₄O₇S

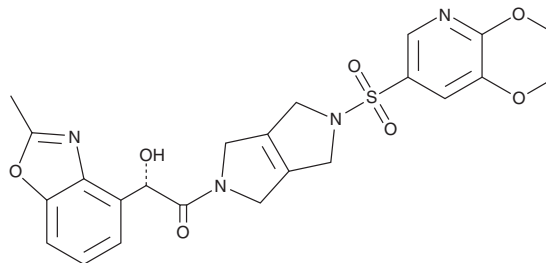
FW: 498.5

Purity: ≥98%

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

FT709 is supplied as a solid. A stock solution may be made by dissolving the FT709 in the solvent of choice, which should be purged with an inert gas. FT709 is soluble in methanol and DMSO.

Description

FT709 is an inhibitor of ubiquitin-specific protease 9X (USP9X; IC₅₀ = 82 nM).¹ It is selective for USP9X over a panel of 20 deubiquitinases (IC₅₀s = >25 μM for all). FT709 (10 μM) decreases the levels of the USP9X substrate centrosomal protein of 55 kDa (CEP55), as well as zinc-finger protein 598 (ZNF598), makorin 1 (MKRN1), and MKRN2, which are involved in the ribosomal quality control pathway, and inhibits the ribosomal stalling response to polyA tracts, a translation quality-control process, in a reporter assay using HEK293 cells.

Reference

1. Clancy, A., Heride, C., Pinto-Fernández, A., *et al.* The deubiquitylase USP9X controls ribosomal stalling. *J. Cell. Biol.* **220**(3), e202004211 (2021).

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, 03/07/2024

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