

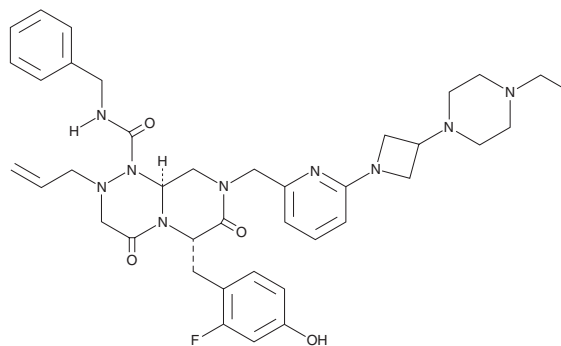
PRODUCT INFORMATION



E-7386

Item No. 46299

CAS Registry No.: 1799824-08-0
Formal Name: (6S,9aS)-8-[[6-[3-(4-ethyl-1-piperazinyl)-1-azetidiny]-2-pyridinyl]methyl]-6-[[2-(2-fluoro-4-hydroxyphenyl)methyl]hexahydro-4,7-dioxo-N-(phenylmethyl)-2-(2-propen-1-yl)-2H-pyrazino[2,1-c][1,2,4]triazine-1(6H)-carboxamide
MF: C₃₉H₄₈FN₉O₄
FW: 725.9
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

E-7386 is supplied as a solid. A stock solution may be made by dissolving the E-7386 in the solvent of choice, which should be purged with an inert gas. E-7386 is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

E-7386 is an inhibitor of the protein-protein interaction between β -catenin and CREB-binding protein (CBP).¹ It inhibits reporter gene expression induced by β -catenin in ECC10 gastric small cell carcinoma cells that express mutant adenomatous polyposis coli and in HEK293 cells (APC; IC₅₀s = 14.7 and 48.4 nM, respectively). E-7386 inhibits the growth of ECC10 cells (IC₅₀ = 4.9 nM) and inhibits tumor growth in an ECC10 mouse xenograft model when administered at a dose of 25-50 mg/kg. It decreases the number of small intestine polyps in an *Apc*^{+/Min} spontaneous colorectal adenoma mouse model and increases intratumoral CD8⁺ T cells and reduces tumor growth in a MMTV-Wnt1 isogenic mouse mammary tumor model.

Reference

1. Yamada, K., Hori, Y., Inoue, S., *et al.* E7386, a selective inhibitor of the interaction between β -Catenin and CBP, Exerts antitumor activity in tumor models with activated canonical Wnt signaling. *Cancer Res.* **81**(4), 1052-1062 (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, 08/05/2026

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