

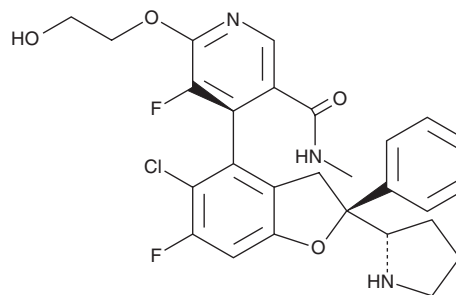
PRODUCT INFORMATION



LAG933

Item No. 45726

CAS Registry No.: 2714434-21-4
Formal Name: 4S-[(2S)-5-chloro-6-fluoro-2,3-dihydro-2-phenyl-2-(2S)-2-pyrrolidinyl-4-benzofuranyl]-5-fluoro-6-(2-hydroxyethoxy)-N-methyl-3-pyridinecarboxamide
MF: C₂₇H₂₆ClF₂N₃O₄
FW: 530.0
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

LAG933 is supplied as a solid. A stock solution may be made by dissolving the LAG933 in the solvent of choice, which should be purged with an inert gas. LAG933 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

LAG933 is an inhibitor of the protein-protein interaction between YAP/TAZ and TEAD (IC₅₀ = 8.6 nM in a TR-FRET assay), downstream effectors of Hippo/MST1/2 signaling, which regulates cell proliferation and is frequently dysregulated in cancer.¹ It reduces growth of Hippo-dependent NCI-H2052 mesothelioma cells and YAP-amplified SF-268 glioma cells (GI₅₀s = 0.04 and 1.11 nM, respectively) but not Hippo-insensitive MKN45 gastric cancer cells (GI₅₀ = >10 nM). Oral administration of LAG933 (30 mg/kg) induces tumor regression in a Hippo-driven MSTO-211H mesothelioma rat xenograft model.

Reference

1. Chapeau, E.A., Sansregret, L., Galli, G.G., *et al.* Direct and selective pharmacological disruption of the YAP-TEAD interface by LAG933 inhibits Hippo-dependent and RAS-MAPK-altered cancers. *Nat. Cancer* **5**(7), 1102-1120 (2024).

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, 06/01/2026

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