

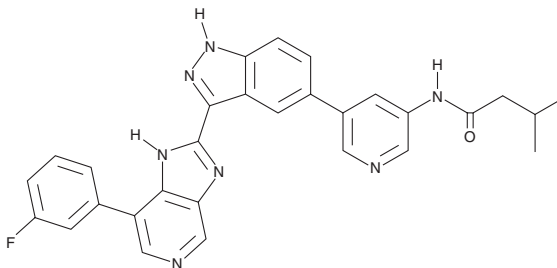
PRODUCT INFORMATION



Adavivint

Item No. 43927

CAS Registry No.: 1467093-03-3
Formal Name: N-[5-[3-[7-(3-fluorophenyl)-3H-imidazo[4,5-c]pyridin-2-yl]-1H-indazol-5-yl]-3-pyridinyl]-3-methyl-butanamide
Synonyms: Lorecivivint, SM-04690
MF: C₂₉H₂₄FN₇O
FW: 505.6
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

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

Description

Adavivint is an inhibitor of Wnt signaling ($EC_{50} = 19.5$ nM in a reporter assay using SW480 cells).¹ It inhibits CMGC family kinases, including CDC-like kinase 2 (Clk2) and dual-specificity tyrosine phosphorylation-regulated kinase 1A (DYRK1A; IC_{50} s = 7.8 and 26.9 nM, respectively), as well as Clk1, Clk3, Clk4, DYRK1B, glycogen synthase kinase 3 β (GSK3 β), and homeodomain-interacting protein kinase 1 (HIPK1), HIPK2, and HIPK3 (IC_{50} s = 16.8-239 nM).² Adavivint also induces activation of YES-associated transcriptional regulator (YAP) in a reporter assay using HEK293A cells ($EC_{50} = 1.2$ nM) in a Clk2-dependent manner.³ It induces chondrogenesis in primary human bone marrow-derived mesenchymal stem cells (MSCs).¹ Adavivint inhibits the replication of pseudorabies virus in infected HeLa, BHK-21, Vero, and PK(15) cells (IC_{50} s = 0.16, 0.19, 0.22, and 0.35 μ M, respectively).⁴ Intra-arterial administration of adavivint (0.3 μ g/animal) decreases disease severity in a surgically induced rat model of osteoarthritis.¹ Adavivint (50 mg/kg) also inhibits tumor growth in a patient-derived organoid xenograft mouse model of colorectal cancer.⁵

References

1. Deshmukh, V., Hu, H., Barroga, C., et al. A small-molecule inhibitor of the Wnt pathway (SM04690) as a potential disease modifying agent for the treatment of osteoarthritis of the knee. *Osteoarthritis Cartilage* **26**(1), 18-27 (2018).
2. Deshmukh, V., O'Green, A.L., Bossard, C., et al. Modulation of the Wnt pathway through inhibition of CLK2 and DYRK1A by lorecivivint as a novel, potentially disease-modifying approach for knee osteoarthritis treatment. *Osteoarthritis Cartilage* **27**(9), 1347-1360 (2019).
3. Bulos, M.L., Grzelak, E.M., Li-Ma, C., et al. Pharmacological inhibition of CLK2 activates YAP by promoting alternative splicing of AMOTL2. *bioRxiv [Preprint]* 1-17 (2023).
4. Wang, C., Wang, T., He, Q., et al. Inhibition of the canonical Wnt/ β -catenin pathway interferes with macropinocytosis to suppress pseudorabies virus proliferation. *Vet. Microbiol.* **301**, 110373 (2025).
5. Xiang, Z., Wang, Y., Ma, X., et al. Targeting the NOTCH2/ADAM10/TCF7L2 axis-mediated transcriptional regulation of Wnt pathway suppresses tumor growth and enhances chemosensitivity in colorectal cancer. *Adv. Sci. (Weinh)* **12**(3), e2405758 (2025).

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, 08/12/2025

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