

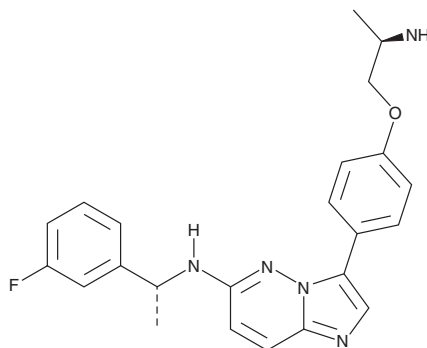
PRODUCT INFORMATION



Taletrectinib

Item No. 45040

CAS Registry No.: 1505514-27-1
Formal Name: 3-[4-[(2R)-2-aminopropoxy] phenyl]-N-[(1R)-1-(3-fluorophenyl)ethyl]-imidazo[1,2-b]pyridazin-6-amine
Synonyms: AB-106, DS-6051b
MF: C₂₃H₂₄FN₅O
FW: 405.5
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

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

Description

Taletrectinib is an inhibitor of c-Ros oncogene 1 (ROS1), tropomyosin-related kinase A (TrkA), and TrkC (IC₅₀s = 0.207, 0.622, and 0.98 nM, respectively).¹ It is selective for these kinases over TrkB (IC₅₀ = 2.28 nM). Taletrectinib (1-10,000 nM) reduces the viability of JFCR-165, JFCR-168, and MGH193-1B patient-derived CD74-ROS1 fusion protein-positive non-small cell lung cancer (NSCLC) cells. *In vivo*, taletrectinib (25-100 mg/kg) reduces tumor volume in Ba/F3 mouse xenograft models expressing CD74-ROS1 wild-type fusion proteins or crizotinib-resistant CD74-ROS1^{G2032R} mutant fusion proteins.

Reference

1. Katayama, R., Gong, B., Togashi, N., *et al.* The new-generation selective ROS1/NTRK inhibitor DS-6051b overcomes crizotinib resistant ROS1-G2032R mutation in preclinical models. *Nat. Commun.* **10**(1), 3604 (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, 02/05/2026

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