

# PRODUCT INFORMATION

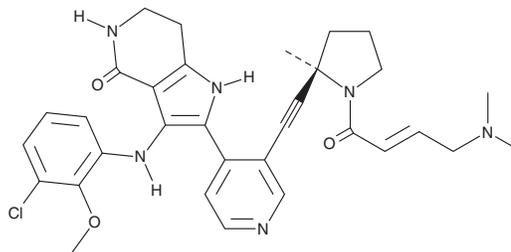


## STX-721

Item No. 44459

**CAS Registry No.:** 2765525-82-2  
**Formal Name:** 3-[(3-chloro-2-methoxyphenyl)amino]-2-[3-[2-[(2R)-1-[(2E)-4-(dimethylamino)-1-oxo-2-buten-1-yl]-2-methyl-2-pyrrolidinyl]ethynyl]-4-pyridinyl]-1,5,6,7-tetrahydro-4H-pyrrolo[3,2-c]pyridin-4-one

**MF:** C<sub>32</sub>H<sub>35</sub>ClN<sub>6</sub>O<sub>3</sub>  
**FW:** 587.1  
**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

STX-721 is supplied as a solid. A stock solution may be made by dissolving the STX-721 in the solvent of choice, which should be purged with an inert gas. STX-721 is slightly soluble (0.1-1 mg/ml) in acetonitrile:water (1:1).

### Description

STX-721 is an irreversible inhibitor of EGFR with an exon 20 insertion (ex20ins) mutation.<sup>1</sup> It selectively inhibits the proliferation of Ba/F3 cells expressing EGFR exon20 A769\_V770 insASV and EGFR exon20 D770\_N771 insSVD (IC<sub>50</sub>s = 5.4 and 5.8 nM, respectively) over those expressing wild-type EGFR (IC<sub>50</sub> = 127.365 nM). STX-721 (50 mg/kg once per day) induces tumor regression in a mouse xenograft model using NCI H2073 non-small cell lung cancer (NSCLC) cells expressing EGFR exon20 D770\_N771 insSVD.

### Reference

1. Milgram, B.C., Borrelli, D.R., Brooijmans, N., *et al.* Discovery of STX-721, a covalent, potent, and highly mutant-selective EGFR/HER2 exon20 insertion inhibitor for the treatment of non-small cell lung cancer. *J. Med. Chem.* **68**(3), 2403-2421 (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, 11/18/2025

#### 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