

PRODUCT INFORMATION

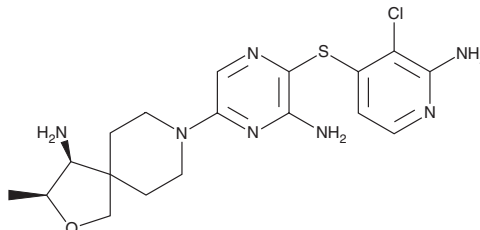


TNO155

Item No. 32525

CAS Registry No.: 1801765-04-7
Formal Name: (3S,4S)-8-[6-amino-5-[(2-amino-3-chloro-4-pyridinyl)thio]-2-pyrazinyl]-3-methyl-2-oxa-8-azaspiro[4.5]decan-4-amine

MF: C₁₈H₂₄ClN₇OS
FW: 421.9
Purity: ≥98%
UV/Vis.: λ_{max}: 224, 269, 355 nm
Supplied as: A crystalline 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

TNO155 is supplied as a crystalline solid. A stock solution may be made by dissolving the TNO155 in the solvent of choice, which should be purged with an inert gas. TNO155 is soluble in DMSO.

Description

TNO155 is an orally bioavailable allosteric inhibitor of Src homology region 2 domain-containing phosphatase 2 (SHP-2; IC₅₀ = 0.011 μM).¹ It is selective for SHP-2 over a panel of 52 kinases and phosphatases (IC₅₀s = >10 μM for all). TNO155 decreases phosphorylation of ERK in, and inhibits proliferation of, KYSE-520 human esophageal cancer cells (IC₅₀s = 0.011 and 0.1 μM, respectively). It acts synergistically with dabrafenib (Item No. 16989) and trametinib (Item No. 16292) to inhibit proliferation of B-RAF^{V600E}-expressing HT-29, RKO, and MDST8 colorectal cancer cells.² TNO155 (2.5, 10, and 20 mg/kg twice per day) reduces tumor growth in a KYSE-520 mouse xenograft model.¹

References

1. LaMarche, M.J., Acker, M., Argintaru, A., *et al.* Identification of TNO155, an allosteric SHP2 inhibitor for the treatment of cancer. *J. Med. Chem.* **63**(22), 13578-13594 (2020).
2. Liu, C., Lu, H., Wang, H., *et al.* Combinations with allosteric SHP2 inhibitor TNO155 to block receptor tyrosine kinase signaling. *Clin. Cancer Res.* **27**(1), 342-354 (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, 10/18/2022

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