

PRODUCT INFORMATION

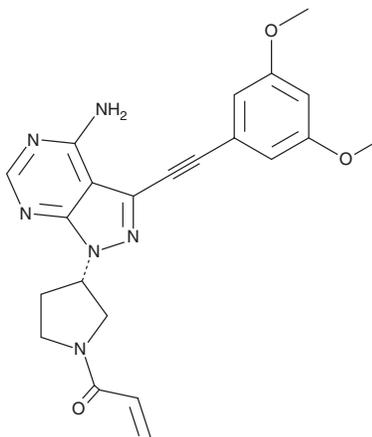


TAS 120

Item No. 21136

CAS Registry No.: 1448169-71-8
Formal Name: 1-[(3S)-3-[4-amino-3-[2-(3,5-dimethoxyphenyl)ethynyl]-1H-pyrazolo[3,4-d]pyrimidin-1-yl]-1-pyrrolidinyl]-2-propen-1-one

MF: C₂₂H₂₂N₆O₃
FW: 418.5
Purity: ≥98%
UV/Vis.: λ_{max}: 210, 301 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

TAS 120 is supplied as a crystalline solid. A stock solution may be made by dissolving the TAS 120 in the solvent of choice, which should be purged with an inert gas. TAS 120 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of TAS 120 in ethanol and DMSO is approximately 1 mg/ml and approximately 30 mg/ml in DMF.

TAS 120 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, TAS 120 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. TAS 120 has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

TAS 120 is an orally bioavailable inhibitor of fibroblast growth factor receptors (FGFRs) that irreversibly inhibits all four FGFR subtypes at nanomolar concentrations.¹ It prevents phosphorylation of FGFRs and impedes downstream signaling pathways in cells.¹ TAS 120 blocks the proliferation of human cancer cell lines with FGFR gene abnormalities but does not alter the growth of cell lines with wild-type FGFR.¹ It inhibits FGFR phosphorylation in tumors and tumor growth in human xenograft mouse models in a dose-dependent manner.¹

Reference

1. Ochiwa, H. Fujita, H., Itoh, K., *et al.* TAS-120, a highly potent and selective irreversible FGFR inhibitor, is effective in tumors harboring various FGFR gene abnormalities. *Mol. Cancer Ther.* **12(11)** Abstract A270, (2013).

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, 12/16/2022

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