

PRODUCT INFORMATION

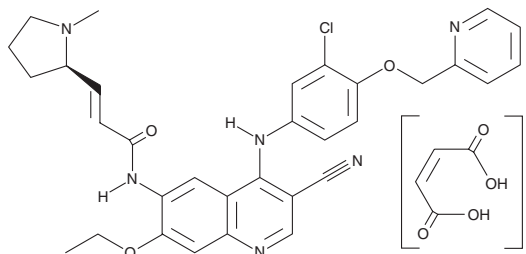


Pyrotinib (maleate)

Item No. 42108

CAS Registry No.: 1397922-61-0
Formal Name: (2E)-N-[4-[[3-chloro-4-(2-pyridinylmethoxy)phenyl]amino]-3-cyano-7-ethoxy-6-quinolinyl]-3-[[[2R]-1-methyl-2-pyrrolidinyl]-2-propenamide, (2Z)-2-butenedioate

Synonyms: SHR-1258
MF: $C_{32}H_{31}ClN_6O_3 \cdot 2C_4H_4O_4$
FW: 815.2
Purity: $\geq 98\%$
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

Pyrotinib (maleate) is supplied as a solid. A stock solution may be made by dissolving the pyrotinib (maleate) in the solvent of choice, which should be purged with an inert gas. Pyrotinib (maleate) is slightly soluble (0.1-1 mg/ml) in acetonitrile and sparingly (1-10 mg/ml) soluble in DMSO.

Pyrotinib (maleate) is slightly soluble (0.1-1 mg/ml) in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Pyrotinib is an inhibitor of EGFR and HER2 (IC_{50} s = 13 and 38 nM, respectively).¹ It selectively inhibits the proliferation of HER2-positive BT474 breast and SKOV3 ovarian cancer cells (IC_{50} s = 5.1 and 43 nM, respectively) over HER2-negative MDA-MB-231 breast cancer cells (IC_{50} = 3,500 nM). *In vivo*, pyrotinib (10 mg/kg per day) decreases tumor growth without affecting body weight in BT474 and SKOV3 mouse xenograft models.

Reference

1. Li, X., Yang, C., Wan, H., *et al.* Discovery and development of pyrotinib: A novel irreversible EGFR/HER2 dual tyrosine kinase inhibitor with favorable safety profiles for the treatment of breast cancer. *Eur. J. Pharm. Sci.* **110**, 51-61 (2017).

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, 10/22/2024

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