

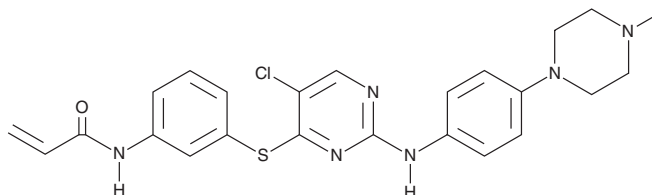
# PRODUCT INFORMATION



**WZ8040**

Item No. 30062

**CAS Registry No.:** 1214265-57-2  
**Formal Name:** N-[3-[[5-chloro-2-[[4-(4-methyl-1-piperazinyl)phenyl]amino]-4-pyrimidinyl]thio]phenyl]-2-propenamide  
**MF:** C<sub>24</sub>H<sub>25</sub>ClN<sub>6</sub>OS  
**FW:** 481.0  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 281 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

WZ8040 is supplied as a crystalline solid. A stock solution may be made by dissolving the WZ8040 in the solvent of choice, which should be purged with an inert gas. WZ8040 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of WZ8040 in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

WZ8040 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, WZ8040 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. WZ8040 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

WZ8040 is an inhibitor of mutant EGFR receptors (EGFRs) with IC<sub>50</sub> values ranging from 2 to 306 nM in Ba/F3 cells.<sup>1</sup> It is selective for EGFR mutants, including the EGFR<sup>Del E746\_A750</sup> and EGFR<sup>L858R</sup> mutations conferring gefitinib sensitivity and the EGFR<sup>Del E746\_A750/T790M</sup> mutation conferring gefitinib resistance (IC<sub>50</sub>s = 2, 6, and 6 nM in Ba/F3 cells), over wild-type EGFR (IC<sub>50</sub> = 1,820 nM in HN11 cells) but also inhibits ERBB2<sup>Ins G776V,C</sup> and wild-type ERBB2 (IC<sub>50</sub>s = 20 and 32 nM, respectively, in Ba/F3 cells). WZ8040 inhibits proliferation of PC-9 cells harboring the EGFR<sup>Del E746\_A750</sup> mutation and gefitinib-resistant PC-9 GR cells harboring the EGFR<sup>T790M</sup> resistant allele (EC<sub>50</sub>s = 6 and 8 nM, respectively).<sup>2</sup>

## References

1. Zhou, W., Ercan, D., Chen, L., *et al.* Novel mutant-selective EGFR kinase inhibitors against EGFR T790M. *Nature* **462**(7276), 1070-1074 (2009).
2. Zhou, W., Ercan, D., Jänne, P.A., *et al.* Discovery of selective irreversible inhibitors for EGFR-T790M. *Bioorg. Med. Chem. Lett.* **21**(2), 638-643 (2011).

### 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, 12/12/2022

## 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](http://WWW.CAYMANCHEM.COM)