

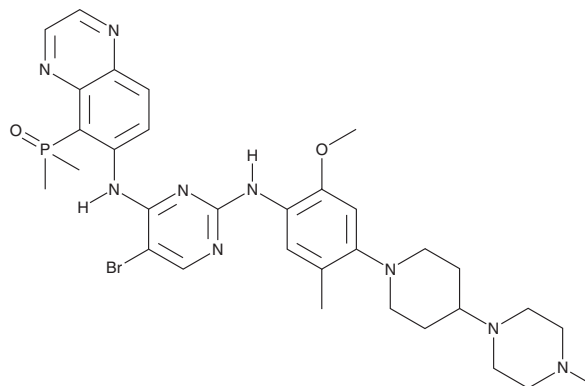
# PRODUCT INFORMATION



## EGFR-IN-7

Item No. 42214

**CAS Registry No.:** 2267329-76-8  
**Formal Name:** 5-bromo-N4-[5-(dimethylphosphinyl)-6-quinoxaliny]-N2-[2-methoxy-5-methyl-4-[4-(4-methyl-1-piperazinyl)-1-piperidinyl]phenyl]-2,4-pyrimidinediamine  
**Synonyms:** EGFR Inhibitor 7, Epidermal Growth Factor Receptor-IN-7, Epidermal Growth Factor Receptor Inhibitor 7, TQB3804  
**MF:**  $C_{32}H_{41}BrN_9O_2P$   
**FW:** 694.6  
**Purity:**  $\geq 95\%$   
**Supplied as:** A solid  
**Storage:**  $-20^{\circ}C$   
**Stability:**  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

EGFR-IN-7 is supplied as a solid. A stock solution may be made by dissolving the EGFR-IN-7 in the solvent of choice, which should be purged with an inert gas. EGFR-IN-7 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

### Description

EGFR-IN-7 is an inhibitor of wild-type and mutant EGFR ( $IC_{50}s = 1.07, 0.46, 0.13, 0.26$ , and  $0.19$  nM for wild-type EGFR, EGFR<sup>d746-750/T790M/C797S</sup>, EGFR<sup>L858R/T790M/C797S</sup>, EGFR<sup>d746-750/T790M</sup>, and EGFR<sup>L858R/T790M</sup>, respectively).<sup>1</sup> It inhibits the proliferation of wild-type EGFR-expressing A431, EGFR<sup>d746-750/T790M/C797S</sup>-expressing Ba/F3, EGFR<sup>d746-750/T790M/C797S</sup>-expressing H1975, and EGFR<sup>d746-750</sup>-expressing PC-9 cells ( $IC_{50}s = 147, 26.8, 163$ , and  $45$  nM, respectively). EGFR-IN-7 reduces tumor growth in triple mutant EGFR mouse xenograft models.

### Reference

1. Liu, X., Zhang, X., Yang, L., *et al.* Abstract 1320: Preclinical evaluation of TQB3804, a potent EGFR C797S inhibitor. *Cancer Res.* 79(13\_Suppl), (2019).

#### 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, 12/18/2024

#### 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