

# PRODUCT INFORMATION



**TX1-85-1**

Item No. 19164

**CAS Registry No.:** 1603845-32-4

**Formal Name:** N-[5-[1-[4-(4-acetyl-1-piperazinyl)cyclohexyl]-4-amino-1H-pyrazolo[3,4-d]pyrimidin-3-yl]-2-phenoxyphenyl]-2-propenamide

**MF:** C<sub>32</sub>H<sub>36</sub>N<sub>8</sub>O<sub>3</sub>

**FW:** 580.7

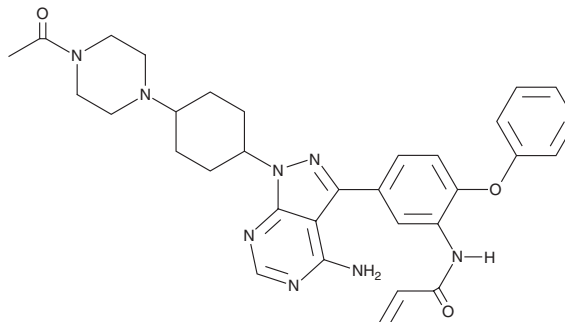
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 266 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

TX1-85-1 is supplied as a crystalline solid. A stock solution may be made by dissolving the TX1-85-1 in the solvent of choice. TX1-85-1 is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of TX1-85-1 in these solvents is approximately 0.2 mg/ml.

## Description

TX1-85-1 is an inhibitor of the epidermal growth factor receptor tyrosine kinase ErbB3 (IC<sub>50</sub> = 23 nM).<sup>1</sup> It also binds to ErbB2, Lyn, and several other Src family kinases in a live cell chemical proteomics assay when used at a concentration of 2 μM. TX1-85-1 inhibits phosphorylation of Akt and ERK1/2 in ErbB3-dependent HCC2935 and OVCAR-8 cells when used at concentrations of 5 and 2 μM, respectively. However, it does not inhibit proliferation of ErbB3-dependent PC9 GR4, HCC827 GR8, or OVCAR-8 cells (EC<sub>50</sub>s = ≥10 μM).

## Reference

1. Xie, T., Lim, S.M., Westover, K.D., *et al.* Pharmacological targeting of the pseudokinase Her3. *Nat. Chem. Biol.* **10**(12), 1006-1012 (2014).

### 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, 11/29/2022

## 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