

# PRODUCT INFORMATION

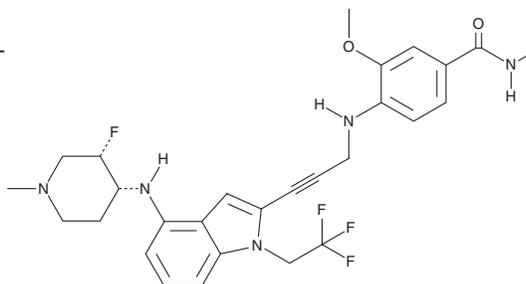


## Rezatapopt

Item No. 44358

**CAS Registry No.:** 2636846-41-6  
**Formal Name:** 4-[[3-[4-[[[(3S,4R)-3-fluoro-1-methyl-4-piperidiny] amino]-1-(2,2,2-trifluoroethyl)-1H-indol-2-yl]-2-propyn-1-yl]amino]-3-methoxy-N-methyl-benzamide

**Synonym:** PC14586  
**MF:** C<sub>28</sub>H<sub>31</sub>F<sub>4</sub>N<sub>5</sub>O<sub>2</sub>  
**FW:** 545.6  
**Purity:** ≥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

Rezatapopt is supplied as a solid. A stock solution may be made by dissolving the rezatapopt in the solvent of choice, which should be purged with an inert gas. Rezatapopt is soluble ( $\geq 10$  mg/ml) in DMSO. Rezatapopt is sparingly soluble (1-10 mg/ml) in ethanol.

### Description

Rezatapopt is a reactivator of p53 containing a tyrosine-to-cysteine substitution at position 220 (p53<sup>Y220C</sup>).<sup>1</sup> It increases p53<sup>Y220C</sup> DNA binding by 1.5-fold in a time-resolved FRET (TR-FRET) assay when used at a concentration of 9 nM. Rezatapopt decreases the proliferation of NUGC-3 gastric carcinoma cells (IC<sub>50</sub> = 504 nM).

### Reference

1. Vu, B.T., Dominique, R., Fahr, B.J., *et al.* Discovery of rezatapopt (PC14586), a first-in-class, small-molecule reactivator of p53 Y220C mutant in development. *ACS Med. Chem. Lett.* **16(1)**, 34-39 (2024).

#### 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, 10/07/2025

#### 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