

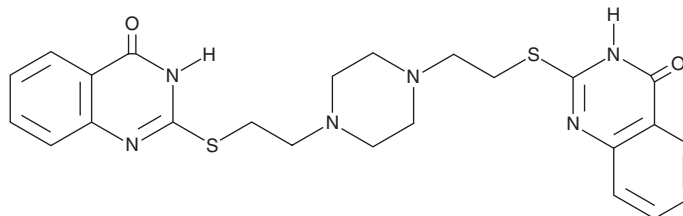
# PRODUCT INFORMATION



**CAY10753**

Item No. 30277

**Formal Name:** 2,2'-((piperazine-1,4-diylbis(ethane-2,1-diyl))bis(sulfanediy))bis(quinazolin-4(3H)-one)  
**MF:** C<sub>24</sub>H<sub>26</sub>N<sub>6</sub>O<sub>2</sub>S<sub>2</sub>  
**FW:** 494.6  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 230, 273 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

CAY10753 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10753 in the solvent of choice, which should be purged with an inert gas. CAY10753 is slightly soluble in DMSO and methanol.

## Description

CAY10753 is an inhibitor of tankyrase 2 (TNKS2; IC<sub>50</sub> = 0.3 nM).<sup>1</sup> It is selective for TNKS2 over TNKS1 (IC<sub>50</sub> = 6.1 nM), as well as poly(ADP-ribose) polymerase 1 (PARP1) and PARP2 (IC<sub>50</sub>s = 89.6 and 37.8 nM, respectively). CAY10753 inhibits proliferation of DLD-1 colorectal cancer cells when used at a concentration of 1 μM.

## Reference

1. Tomassi, S., Pfahler, J., Mautone, N., *et al.* From PARP1 to TNKS2 inhibition: A structure-based approach. *Med. Chem. Lett.* (2020).

### 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, 11/07/2025

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