

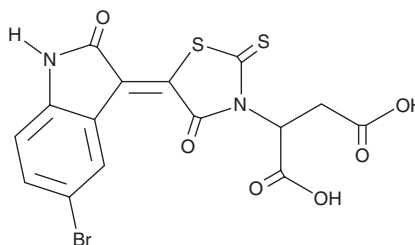
# PRODUCT INFORMATION



**CID-5721353**

Item No. 22967

**CAS Registry No.:** 301356-95-6  
**Formal Name:** 2-[5-(5-bromo-1,2-dihydro-2-oxo-3H-indol-3-ylidene)-4-oxo-2-thioxo-3-thiazolidinyl]-butanedioic acid  
**MF:** C<sub>15</sub>H<sub>9</sub>BrN<sub>2</sub>O<sub>6</sub>S<sub>2</sub>  
**FW:** 457.3  
**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

CID-5721353 is supplied as a solid. A stock solution may be made by dissolving the CID-5721353 in the solvent of choice, which should be purged with an inert gas. CID-5721353 is soluble in organic solvents such as acetonitrile and DMSO.

## Description

CID-5721353 is an inhibitor of the transcriptional repressor B cell lymphoma 6 (Bcl-6;  $K_i = 147 \mu\text{M}$ ).<sup>1</sup> It is selective for Bcl-6 over Kaiso, hypermethylated in cancer 1 (HIC1), and promyelocytic leukemia zinc finger (PLZF) at 50  $\mu\text{M}$ . CID-5721353 is cytotoxic against Bcl-6-dependent SU-DLH-6, SU-DLH-4, Farage, OCI-LY7, OCI-LY1, and OCI-LY10 diffuse large B cell lymphoma (DLBCL) cells ( $GI_{50} = 0.024\text{-}0.936 \text{ mM}$ ) but not Bcl-6-independent Toledo or OCI-LY4 DLBCL cells ( $GI_{50} = >15 \text{ mM}$  for both). It increases expression of the Bcl-6 target genes *ATR*, *TP53*, *CD69*, *p21*, and *CD44* in SU-DLH-6 and SU-DHL-4 cells when used at a concentration of 50  $\mu\text{M}$ . CID-5721353 (50 mg/kg) reduces tumor growth in OCI-LY7 and SU-DHL-6 mouse xenograft models.

## Reference

1. Cerchietti, L.C., Ghetu, A.F., Zhu, X., *et al.* A small-molecule inhibitor of BCL6 kills DLBCL cells in vitro and in vivo. *Cancer Cell* **17**(4), 400-411 (2010).

**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, 01/18/2024

## 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