

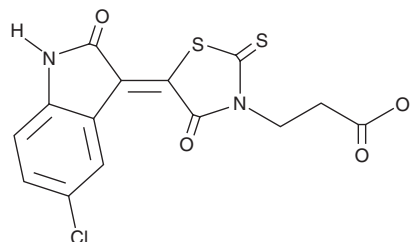
# PRODUCT INFORMATION



## FX1

Item No. 23908

**CAS Registry No.:** 1426138-42-2  
**Formal Name:** (5Z)-5-(5-chloro-1,2-dihydro-2-oxo-3H-indol-3-ylidene)-4-oxo-2-thioxo-3-thiazolidinepropanoic acid  
**MF:** C<sub>14</sub>H<sub>9</sub>ClN<sub>2</sub>O<sub>4</sub>S<sub>2</sub>  
**FW:** 368.8  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 232, 276, 416 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

FX1 is supplied as a crystalline solid. A stock solution may be made by dissolving the FX1 in the solvent of choice. FX1 is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of FX1 in these solvents is approximately 30 mg/ml.

### Description

FX1 is an inhibitor of the transcription factor B cell lymphoma 6 (BCL6; IC<sub>50</sub> = ~35 μM).<sup>1</sup> It is selective for BCL6 over a panel of 50 kinases at a concentration of 10 μM. FX1 (50 μM) reduces recruitment of the BCL6 corepressor proteins SMRT and BCOR to the BCL6 target loci *CDKN1A*, *CD69*, and *CXCR4* in a ChIP assay. It represses growth of Farage, SU-DLH-6, DOHH-2, OCI-Ly7, RCK8, SU-DHL-4, OCI-Ly1, and SC-1 BCL6-dependent cell lines (GI<sub>50</sub>s = 16-54 μM) but not OCI-LY1-B50, Karpas 433, OCI-Ly19, or Toledo BCL6-independent cell lines (GI<sub>50</sub>s = >125 μM). *In vivo*, FX1 (50 mg/kg) reduces tumor volume in an HBL-1 large B cell lymphoma mouse xenograft model.

### Reference

1. Cardenas, M.G., Yu, W., Beguelin, W., *et al.* Rationally designed BCL6 inhibitors target activated B cell diffuse large B cell lymphoma. *J. Clin. Invest.* **126**(9), 3351-3362 (2016).

#### 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, 12/22/2022

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