

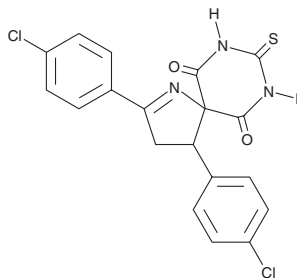
# PRODUCT INFORMATION



## SCR130

Item No. 37779

**CAS Registry No.:** 2377858-38-1  
**Formal Name:** 2,4-bis(4-chlorophenyl)-8-thioxo-1,7,9-triazaspiro[4.5]dec-1-ene-6,10-dione  
**MF:** C<sub>19</sub>H<sub>13</sub>Cl<sub>2</sub>N<sub>3</sub>O<sub>2</sub>S  
**FW:** 418.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 249, 292 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

SCR130 is supplied as a crystalline solid. A stock solution may be made by dissolving the SCR130 in the solvent of choice, which should be purged with an inert gas. SCR130 is soluble in DMSO (warmed, sonicated).

### Description

SCR130 is an inhibitor of non-homologous end joining (NHEJ) and a derivative of the DNA ligase IV inhibitor SCR7 (Item No. 18015).<sup>1</sup> It selectively inhibits DNA ligase IV- over ligase I- or ligase III-mediated end-joining in cell-free assays when used at a concentration of 200 μM. SCR130 is cytotoxic to NALM6 and Reh leukemia and HeLa cervical cancer cells (IC<sub>50</sub>s = 2.2, 14.1, and 5.9 μM, respectively) and induces DNA double-strand breaks (DSBs) in Reh cells when used at concentrations ranging from 7 to 21 μM.

### Reference

1. Ray, U., Raul, S.K., Gopinatha, V.K., *et al.* Identification and characterization of novel SCR7-based small-molecule inhibitor of DNA end-joining, SCR130 and its relevance in cancer therapeutics. *Mol. Carcinog.* **59(6)**, 618-628 (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, 07/11/2023

#### 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