

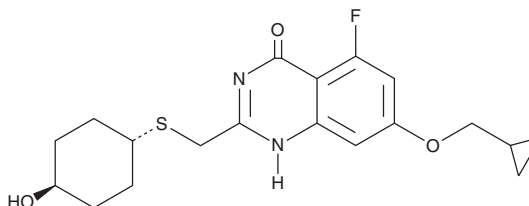
# PRODUCT INFORMATION



**RBN-012759**

Item No. 36137

**CAS Registry No.:** 2360851-29-0  
**Formal Name:** 7-(cyclopropylmethoxy)-5-fluoro-2-[[[(trans-4-hydroxycyclohexyl)thio]methyl]-4(3H)-quinazolinone  
**MF:** C<sub>19</sub>H<sub>23</sub>FN<sub>2</sub>O<sub>3</sub>S  
**FW:** 378.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 243 nm  
**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

RBN-012759 is supplied as a solid. A stock solution may be made by dissolving the RBN-012759 in the solvent of choice, which should be purged with an inert gas. RBN-012759 is soluble in DMSO.

## Description

RBN-012759 is an inhibitor of poly(ADP-ribose) polymerase 14 (PARP14; IC<sub>50</sub> = 0.003 μM).<sup>1</sup> It is selective for PARP14 over PARP1-4, -5a, -5b, -6-12, -15, and -16 (IC<sub>50</sub>s = 1->100 μM). RBN-012759 (10 μM) inhibits IFN-γ-induced PARP14 auto-MARylation in CFPAC-1 cells. It inhibits IL-4-induced protumor M2-like macrophage polarization in isolated human primary macrophages. RBN-012759 (1 μM) also induces expression of the genes encoding the pro-inflammatory chemokines chemokine (C-C motif) ligand 13 (CCL13), CCL19, and chemokine (C-X-C motif) ligand 11 (CXCL11) in human kidney cancer tumor explants.

## Reference

1. Schenkel, L.B., Molina, J.R., Swinger, K.K., *et al.* A potent and selective PARP14 inhibitor decreases protumor macrophage gene expression and elicits inflammatory responses in tumor explants. *Cell Chem. Biol.* **28(8)**, 1158-1168.e13 (2021).

### 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, 11/02/2022

## 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