

PRODUCT INFORMATION

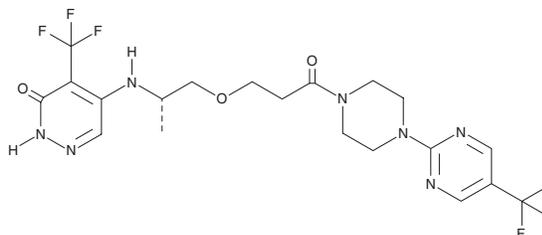


RBN-2397

Item No. 33558

CAS Registry No.: 2381037-82-5
Formal Name: 5-[[[(1S)-1-methyl-2-[3-oxo-3-[4-[5-(trifluoromethyl)-2-pyrimidinyl]-1-piperazinyl]propoxy]ethyl]amino]-4-(trifluoromethyl)-3(2H)-pyridazinone

MF: C₂₀H₂₃F₆N₇O₃
FW: 523.4
Purity: ≥98%
UV/Vis.: λ_{max}: 221, 254 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

RBN-2397 is supplied as a crystalline solid. A stock solution may be made by dissolving the RBN-2397 in the solvent of choice, which should be purged with an inert gas. RBN-2397 is soluble in DMSO and dimethyl formamide. The solubility of RBN-2397 in these solvents is approximately 1 mg/ml, respectively.

Description

RBN-2397 is an inhibitor of poly(ADP-ribose) polymerase 7 (PARP7).¹ It stabilizes wild-type PARP7 levels, but not levels of the catalytically inactive mutant PARP7^{H532A}, in COS-1 cells expressing wild-type or mutant PARP7, respectively, when used at a concentration of 100 nM. RBN-2397 (100 nM) also decreases wild-type PARP7 catalytic activity in the same cells.

Reference

1. Rasmussen, M., Tan, S., Somisetty, V.S., *et al.* PARP7 and mono-ADP-ribosylation negatively regulate estrogen receptor α signaling in human breast cancer cells. *Cells* **10(3)**, 623 (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, 11/08/2022

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