

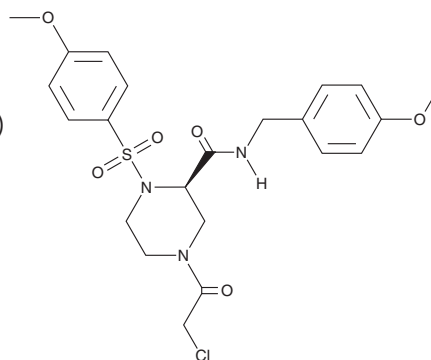
PRODUCT INFORMATION



(R)-SKBG-1

Item No. 42091

CAS Registry No.: 2955634-67-8
Formal Name: (2R)-4-(2-chloroacetyl)-N-[(4-methoxyphenyl)methyl]-1-[(4-methoxyphenyl)sulfonyl]-2-piperazinecarboxamide
MF: C₂₂H₂₆ClN₃O₆S
FW: 496.0
Purity: ≥95%
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

(R)-SKBG-1 is supplied as a solid. A stock solution may be made by dissolving the (R)-SKBG-1 in the solvent of choice, which should be purged with an inert gas. (R)-SKBG-1 is soluble in the organic solvent DMSO.

Description

(R)-SKBG-1 is a degrader of mRNA encoding the androgen receptor (AR).¹ It decreases mRNA levels of AR and AR splice variant seven (AR-V7) and protein levels of AR and AR-V7 in 22Rv1 prostate cancer cells when used at a concentration of 20 μM. (R)-SKBG-1 (20 μM) reduces protein levels of AR and AR-V7 in wild-type 22Rv1 cells but not in *NONO*^{-/-} 22Rv1 cells, which lack the RNA-binding protein non-POU domain-containing octamer-binding protein (NONO). It increases the number of NONO nuclear foci and NONO-bound immature mRNA transcripts in 22Rv1 cells. (R)-SKBG-1 is cytotoxic to 22Rv1 cells (IC₅₀ = 44 μM).

Reference

1. Kathman, S.G., Koo, S.J., Lindsey, G.L., *et al.* Remodeling oncogenic transcriptomes by small molecules targeting NONO. *Nat. Chem. Biol.* **19**(7), 825-836 (2023).

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, 09/11/2024

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