

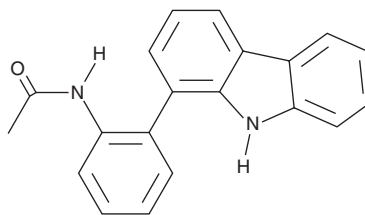
PRODUCT INFORMATION



GeA-69

Item No. 39533

CAS Registry No.: 2143475-98-1
Formal Name: N-[2-(9H-carbazol-1-yl)phenyl]-acetamide
MF: C₂₀H₁₆N₂O
FW: 300.4
Purity: ≥98%
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

GeA-69 is supplied as a solid. A stock solution may be made by dissolving the GeA-69 in the solvent of choice, which should be purged with an inert gas. GeA-69 is soluble in methanol and DMSO.

Description

GeA-69 is an inhibitor of poly(ADP-ribose) polymerase 14 macrodomain 2 (PARP14 MD2).¹ It binds to PARP14 MD2 (K_d = 860 nM) and inhibits PARP14 MD2 relocalization to the sites of DNA damage in U2OS cells when used at a concentration of 50 μM.

Reference

1. Schuller, M., Riedel, K., Gibbs-Seymour, I., *et al.* Discovery of a selective allosteric inhibitor targeting macrodomain 2 of polyadenosine-diphosphate-ribose polymerase 14. *ACS Chem. Biol.* **12**(11), 2866-2874 (2017).

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, 10/26/2023

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