

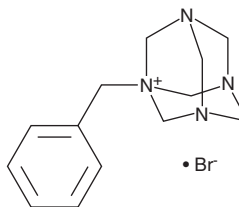
PRODUCT INFORMATION



Roslin-2

Item No. 21985

CAS Registry No.: 29574-21-8
Formal Name: 1-(phenylmethyl)-3,5,7-triaza-1-azoniatricyclo[3.3.1.1^{3,7}]decane, monobromide
MF: C₁₃H₁₉N₄ • Br
FW: 311.2
Purity: ≥98%
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

Roslin-2 is supplied as a crystalline solid. A stock solution may be made by dissolving the roslin-2 in the solvent of choice, which should be purged with an inert gas. Roslin-2 is soluble in the organic solvent DMSO.

Description

Roslin-2 is a reactivator of p53.¹ Roslin-2 (111, 333, and 1000 μM) impairs the interaction between focal adhesion kinase (FAK) and p53 and restores p53 transcriptional activity. It reverses FAK-mediated suppression of the expression of p53-dependent p21, MDM2, and Bax in wild-type p53 HCT116 colon and MCF-7 breast cancer cells, but not in p53 mutant SW620 colon cancer cells, when used at concentrations ranging from 1 to 50 μM. Roslin-2 (10 μM) potentiates the pro-apoptotic effect of doxorubicin (Item No. 15007) and 5-fluorouracil (Item No. 14416) in p53^{+/+}, but not p53^{-/-}, HCT116 cells. It reduces tumor growth and activates p21 in p53^{+/+}, but not p53^{-/-} tumors, in HCT116 mouse xenograft models when administered at a dose of 60 mg/kg.

Reference

1. Golubovskaya, V.M., Ho, B., Zheng, M., *et al.* Disruption of focal adhesion kinase and p53 interaction with small molecule compound R2 reactivated p53 and blocked tumor growth. *BMC Cancer* **13:342**, (2013).

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, 11/14/2022

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