

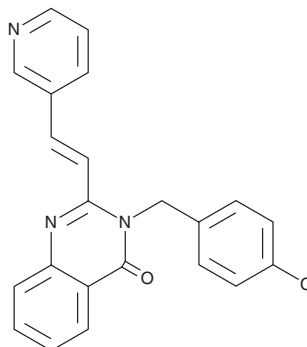
PRODUCT INFORMATION



RAD51-IN-17

Item No. 42253

CAS Registry No.: 2101739-18-6
Formal Name: 3-[(4-chlorophenyl)methyl]-2-[(1E)-2-(3-pyridinyl)ethenyl]-4(3H)-quinazolinone
Synonyms: AFN39186, RAD51 Inhibitor 17
MF: C₂₂H₁₆ClN₃O
FW: 373.8
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

RAD51-IN-17 is supplied as a solid. A stock solution may be made by dissolving the RAD51-IN-17 in the solvent of choice, which should be purged with an inert gas. RAD51-IN-17 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

RAD51-IN-17 is an inhibitor of RAD51 recombinase.¹ It inhibits DNA damage-induced RAD51 foci formation in MDA-MB-231 triple-negative breast cancer (TNBC) cells when used at a concentration of 10 μM. RAD51-IN-17 (10 μM) enhances irradiation-induced reductions in the proliferation of MDA-MB-231 cells.

Reference

1. Ward, A., Dong, L., Harris, J.M., *et al.* Quinazolinone derivatives as inhibitors of homologous recombinase RAD51. *Bioorg. Med. Chem. Lett.* **27**(14), 3096-3100. (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, 11/12/2024

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