

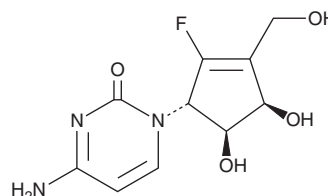
PRODUCT INFORMATION



RX-3117

Item No. 27922

CAS Registry No.: 865838-26-2
Formal Name: 4-amino-1-[(1S,4R,5S)-2-fluoro-4,5-dihydroxy-3-(hydroxymethyl)-2-cyclopenten-1-yl]-2(1H)-pyrimidinone
Synonyms: Fluorocyclopentenylcytosine, TV-1360
MF: C₁₀H₁₂FN₃O₄
FW: 257.2
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

RX-3117 is supplied as a solid. A stock solution may be made by dissolving the RX-3117 in the solvent of choice, which should be purged with an inert gas. RX-3117 is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

RX-3177 is a nucleoside derivative and an anticancer agent.¹ It is cytotoxic to HCT116 colorectal, MDA-MB-231 breast, PANC-1 pancreatic, A549 lung, and U251 brain cancer cells (IC₅₀s = 0.39, 0.18, 0.62, 0.34, and 0.83 μM, respectively). RX-3177 (1 and 10 μM) induces DNA damage and apoptosis in A2780 ovarian cancer cells. *In vivo*, RX-3177 (50-500 mg/kg) reduces mean tumor volume in HCT116, HT-29, Caki-1, MIA PaCa-2, and BxPC-3 mouse xenograft models.²

References

1. Balboni, B., El Hassouni, B., Honeywell, R.J., *et al.* RX-3117 (fluorocyclopentenyl cytosine): A novel specific antimetabolite for selective cancer treatment. *Expert Opin. Investig. Drugs* **28**(4), 311-312 (2019).
2. Yang, M.Y., Lee, Y.B., Ahn, C.-H., *et al.* A novel cytidine analog, RX-3117, shows potent efficacy in xenograft models, even in tumors that are resistant to gemcitabine. *Anticancer Res.* **34**(12), 6951-6959 (2014).

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, 09/25/2024

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