

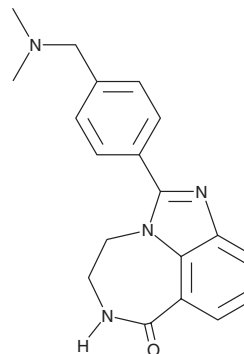
PRODUCT INFORMATION



AG-14361

Item No. 24677

CAS Registry No.: 328543-09-5
Formal Name: 2-[4-[(dimethylamino)methyl]phenyl]-5,6-dihydroimidazo[4,5,1-jk][1,4]benzodiazepin-7(4H)-one
MF: C₁₉H₂₀N₄O
FW: 320.4
Purity: ≥98%
UV/Vis.: λ_{max}: 210, 240, 312 nm
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

AG-14361 is supplied as a crystalline solid. A stock solution may be made by dissolving the AG-14361 in the solvent of choice. AG-14361 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of AG-14361 in these solvents is approximately 2, 20, and 10 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of AG-14361 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of AG-14361 in PBS, pH 7.2, is approximately 0.2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

AG-14361 is an inhibitor of poly(ADP-ribose) polymerase 1 (PARP1; K_i = 6.3 nM for the human enzyme).¹ It inhibits PARP1 activity in permeabilized and intact SW620 cells (IC₅₀s = 29 and 14 nM, respectively).² AG-14361 (0.4 μM) potentiates the growth inhibitory effects of temozolomide (Item No. 14163) in HCT116, A2480, and CP70 cells and of the topoisomerase I inhibitor topotecan (Item No. 14129) by greater than three-fold in PARP1^{+/+} primary embryonic fibroblasts.^{3,4} It slows the rejoining of camptothecin-induced DNA strand breaks.⁴ AG-14361 also inhibits cell growth in a PARP1-independent manner with GI₅₀ values of 66 and 65 μM, respectively, in cell lines derived from PARP1^{+/+} and PARP1^{-/-} mice.² AG-14361 (15 mg/kg), in combination with temozolomide, leads to tumor regression in an SW620 mouse xenograft model.

References

1. Skalitzky, D.J., Marakovits, J.T., Maegley, K.A., *et al. J. Med. Chem.* **46**(2), 210-213 (2003).
2. Calabrese, C.R., Almassy, R., Barton, S., *et al. J. Natl. Cancer Inst.* **96**(1), 56-67 (2004).
3. Curtin, N.J., Wang, L.Z., Yiakouvaki, A., *et al. Clin. Cancer Res.* **10**(3), 881-889 (2004).
4. Smith, L.M., Willmore, E., Austin, C.A., *et al. Clin. Cancer Res.* **11**(23), 8449-8457 (2005).

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, 12/19/2022

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