

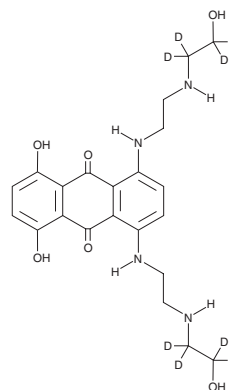
PRODUCT INFORMATION



Mitoxantrone-d₈

Item No. 29414

CAS Registry No.: 1189974-82-0
Formal Name: 1,4-dihydroxy-5,8-bis[[2-[(2-hydroxyethyl-1,1,2,2-d₄)amino]ethyl]amino]-9,10-anthracenedione
MF: C₂₂H₂₀D₈N₄O₆
FW: 452.5
Chemical Purity: ≥95% (Mitoxantrone)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₈); ≤1% d₀
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

Mitoxantrone-d₈ is intended for use as an internal standard for the quantification of mitoxantrone (Item No. 14842) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Mitoxantrone-d₈ is supplied as a solid. A stock solution may be made by dissolving the mitoxantrone-d₈ in the solvent of choice, which should be purged with an inert gas. Mitoxantrone-d₈ is soluble in organic solvents such as methanol and DMSO.

Description

Mitoxantrone is an inhibitor of DNA topoisomerase II α and HIV-1 integrase (IC₅₀s = 5.3 and 3.8 μ M, respectively).^{1,2} It intercalates into DNA in a cell-free assay when used at a concentration of 2 μ M, as well as inhibits DNA synthesis and induces DNA-protein crosslinks *in vitro* in a concentration-dependent manner.^{1,3} Mitoxantrone inhibits the growth of K562 leukemia and K562-derived etoposide-resistant K/VP.5 cells (IC₅₀s = 0.42 and 1.68 μ M, respectively).¹ It also inhibits the membrane fusion of a mixed population of HEK293T cells expressing either angiotensin-converting enzyme 2 (ACE2) or the extracellular domain of severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) spike glycoprotein, also known as surface glycoprotein.⁴ Mitoxantrone (0.2 μ M) inhibits the entry of SARS-CoV-2 into Vero E6 cells. Formulations containing mitoxantrone have been used in the treatment of cancer and multiple sclerosis.

References

1. Hasinoff, B.B., Liang, H., Wu, X., *et al.* The structure-based design, synthesis and biological evaluation of DNA-binding bisintercalating bisanthrapyrazole anticancer compounds. *Bioorg. Med. Chem.* **16**(7), 3959-3968 (2008).
2. Carlson, H.A., Masukawa, K.M., Rubins, K., *et al.* Developing a dynamic pharmacophore model for HIV-1 integrase. *J. Med. Chem.* **43**(111), 2100-2114 (2000).
3. Fox, M.E. and Smith, P.J. Long-term inhibition of DNA synthesis and the persistence of trapped topoisomerase II complexes in determining the toxicity of the antitumor DNA intercalators mAMSA and mitoxantrone. *Cancer Res.* **50**(18), 5813-5818 (1990).
4. Zhang, Q., Radvak, P., Lee, J., *et al.* Mitoxantrone modulates a heparan sulfate-spike complex to inhibit SARS-CoV-2 infection. *Sci. Rep.* **12**(1), 6294 (2022).

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, 12/12/2022

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