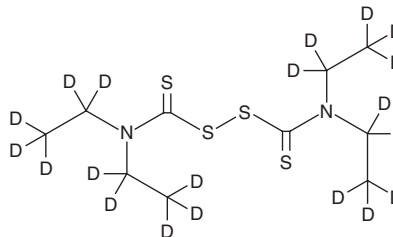


PRODUCT INFORMATION



Disulfiram-d₂₀ Item No. 28536

CAS Registry No.: 1216403-88-1
Formal Name: bis(1,1,2,2,2-pentadeuterioethyl) carbamothioylsulfanyl N,N-bis(1,1,2,2,2-pentadeuterioethyl)carbamodithioate
Synonym: Tetraethylthiuram disulfide-d₂₀
MF: C₁₀D₂₀N₂S₄
FW: 316.7
Chemical Purity: ≥98% (Disulfiram)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₂₀); ≤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

Disulfiram-d₂₀ is intended for use as an internal standard for the quantification of disulfiram (Item No. 15303) 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).

Disulfiram-d₂₀ is supplied as a solid. A stock solution may be made by dissolving the disulfiram-d₂₀ in the solvent of choice, which should be purged with an inert gas. Disulfiram-d₂₀ is soluble in methanol, DMSO, and dimethyl formamide.

Description

Disulfiram is a copper and zinc chelator and an irreversible inhibitor of aldehyde dehydrogenase (IC₅₀ = 0.1 mM).¹ It also inhibits the copper-dependent enzyme dopamine β-hydroxylase, which prevents the breakdown of dopamine.² When in complex with copper, disulfiram has been shown to inhibit purified 20S proteasome (IC₅₀ = 7.5 μM) and 26S proteasome (IC₅₀ = 20 μM) from MDA-MB-0231 breast cancer cells.^{3,4} Disulfiram (at 250 nM) induces reactive oxygen species, activates JNK and p38 pathways, and inhibits NF-κB activity, which suppresses self-renewal in cancer stem cells.^{5,6}

References

1. Kraemer, R.J. and Deitrich, R.A. Isolation and characterization of human liver aldehyde dehydrogenase. *J. Biol Chem.* **243(24)**, 6402-6408 (1968).
2. Haile, C.N., De La Garza, R., II, Mahoney, J.J., III, *et al.* The impact of disulfiram treatment on the reinforcing effects of cocaine: A randomized clinical trial. *PLoS One* **7(11)**, e47702 (2012).
3. Chen, D., Cui, Q.C., Yang, H., *et al.* Disulfiram, a clinically used anti-alcoholism drug and copper-binding agent, induces apoptotic cell death in breast cancer cultures and xenografts via inhibition of the proteasome activity. *Cancer Res.* **66(21)**, 10425-10433 (2006).
4. Schmitt, S.M., Frezza, M., and Dou, Q.P. New applications of old metal-binding drugs in the treatment of human cancer. *Front. Biosci. (Schol. Ed.)* **4**, 375-391 (2012).
5. Liu, P., Brown, S., Goktug, T., *et al.* Cytotoxic effect of disulfiram/copper on human glioblastoma cell lines and ALDH-positive cancer-stem-like cells. *Br. J. Cancer* **107(9)**, 1488-1497 (2012).
6. Chiba, T., Suzuki, E., Yuki, K., *et al.* Disulfiram eradicates tumor-initiating hepatocellular carcinoma cells in ROS-p38 MAPK pathway-dependent and -independent manners. *PLoS One* **9(1)**, e84807 (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, 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