PRODUCT INFORMATION



α-Hydroxyglutaric Acid

Item No. 25894

CAS Registry No.:	2889-31-8	
Formal Name:	2-hydroxy-pentanedioic acid	
Synonyms:	2-HG, 2-Hydroxyglutaric Acid	
MF:	C ₅ H ₈ O ₅	O O
FW:	148.1	но он
Purity:	≥95%	
Supplied as:	A solution in ethanol	ÓH
Storage:	-20°C	
Stability:	≥1 year	

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

 α -Hydroxyglutaric acid (2-HG) is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of 2-HG in these solvents is approximately 30 mg/ml.

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. If an organic solvent-free solution of 2-HG is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of 2-HG in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

2-HG is an α -hydroxy acid. It is metabolized to 2-oxoglutarate by D- and L-2-hydroxyglutarate dehydrogenases, and mutations in these enzymes lead to 2-hydroxyglutaric aciduria, a neurometabolic disorder characterized by increased levels of L-2-hydroxyglutaric acid (Item No. 21123).¹⁻³ 2-HG can also be metabolized from isocitrate by mutated forms of isocitrate dehydrogenase 1 (IDH1) and IDH2.^{4,5} 2-HG is structurally similar to α -ketoglutarate, the product of wild-type isocitrate dehydrogenases, and competitively inhibits α-ketoglutarate-dependent dioxygenases, including histone lysine demethylases and DNA hydroxylases.5-7

References

- 1. Rzem, R., Veiga-da-Cunha, M., Noël, G., et al. Proc. Natl. Acad. Sci. U.S.A. 101(48), 16849-16854 (2004).
- 2. Struys, E.A., Verhoeven, N.M., Roos, B., et al. Clin. Chem. 49(7), 1133-1138 (2003).
- 3. Struys, E.A., Salomons, G.S., Achouri, Y., et al. Am. J. Hum. Genet. 76(2), 358-360 (2005).
- 4. Ward, P., Patel, J., Wise, D.R., et al. Cancer Cell 17(3), 225-234 (2010).
- 5. Yang, H., Ye, D., Guan, K.L., et al. Clin. Cancer Res. 18(20), 5562-5571 (2012).
- 6. Xu, W., Yang, H., Liu, Y., et al. Cancer Cell 19(1), 17-30 (2011).
- 7. Chowdhury, R., Yeoh, K.K., Tian, Y.M., et al. EMBO Rep. 12(5), 463-469 (2011).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFFTY 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

uyer 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, 10/23/2018

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