

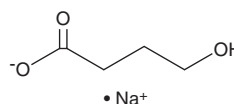
PRODUCT INFORMATION



GHB (sodium salt)

Item No. 9002506

CAS Registry No.: 502-85-2
Formal Name: 4-hydroxy-butanoic acid, monosodium salt
Synonyms: γ -Hydroxybutyric Acid, NSC 84223
MF: $C_4H_7O_3 \cdot Na$
FW: 126.1
Purity: $\geq 95\%$
Supplied as: A crystalline solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



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

Description

GHB (sodium salt) (Item No. 9002506) is an analytical reference standard categorized as a CNS depressant.¹ GHB is a metabolite of γ -butyrolactone (Item No. 18706).² It has a high potential for abuse.^{3,4} GHB is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

References

1. Busardò, F.P. and Kyriakou, C. GHB in biological specimens: Which cut-off levels should be taken into consideration in forensic toxicological investigation? *Recent Pat. Biotechnol.* **8(3)**, 206-214 (2014).
2. Winstock, A.R. and Mitcheson, L. New recreational drugs and the primary care approach to patients who use them. *BMJ* **344**, 1-10 (2012).
3. Carter, L.P., Koek, W., and France, C.P. Behavioral analyses of GHB: Receptor mechanisms. *Pharmacol. Ther.* **121(1)**, 100-114 (2009).
4. Griffiths, R.R. and Johnson, M.W. Relative abuse liability of hypnotic drugs: A conceptual framework and algorithm for differentiating among compounds. *J. Clin. Psychiatry* **66(Suppl 9)**, 31-41 (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, 07/13/2023

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