

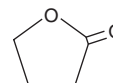
PRODUCT INFORMATION



γ -Butyrolactone

Item No. 18706

CAS Registry No.: 96-48-0
Formal Name: dihydro-2(3H)-furanone
Synonyms: GBL, γ -Hydroxybutyric Acid lactone, NIH 10540, NSC 4592
MF: C₄H₆O₂
FW: 86.1
Purity: \geq 95%
Supplied as: A solution in acetonitrile
Storage: -20°C
Stability: \geq 5 years



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

Description

γ -Butyrolactone (Item No. 18706) is an analytical reference standard categorized as a prodrug form of GHB (Item No. 9002506).^{1,2} γ -Butyrolactone is commonly used as a solvent and reagent in chemistry applications.¹⁻³ γ -Butyrolactone is regulated as a List I chemical in the United States. This product is intended for research and forensic applications.

References

1. Winstock, A.R. and Mitcheson, L. New recreational drugs and the primary care approach to patients who use them. *BMJ* **344**, 1-10 (2012).
2. Busardó, F.P. and Jones, A.W. GHB pharmacology and toxicology: Acute intoxication, concentrations in blood and urine in forensic cases and treatment of the withdrawal syndrome. *Curr. Neuropharmacol.* **13(1)**, 47-70 (2015).
3. U.S. Department of Justice GHB Analogs, National Drug Intelligence Center, in Information Bulletin, 2002-L0424-003, (2002).

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, 06/16/2023

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