

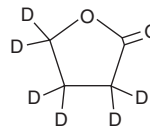
PRODUCT INFORMATION



γ -Butyrolactone-d₆

Item No. 31373

CAS Registry No.: 77568-65-1
Formal Name: dihydro-d₂-2(3H)-furanone-3,3,4,5-d₄
Synonyms: GBL-d₆, γ -Hydroxybutyric Acid lactone-d₆
MF: C₄D₆O₂
FW: 92.1
Chemical Purity: ≥98% (γ -Butyrolactone)
Deuterium
Incorporation: ≥99% deuterated forms (d₁-d₆); ≤1% d₀
UV/Vis.: λ_{max} : 210 nm
Supplied as: A solution in acetonitrile
Storage: -20°C
Stability: ≥4 years



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

Description

γ -Butyrolactone-d₆ (Item No. 31373) is intended for use as an internal standard for the quantification of γ -butyrolactone (Item No. 18706) 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).

γ -Butyrolactone is a prodrug form of GHB (Item No. 9002506).¹ γ -Butyrolactone is commonly used as a solvent and reagent in chemistry applications.¹⁻³ γ -Butyrolactone-d₆ is regulated as a List I chemical in the United States. This product is intended for research and forensic applications.

References

1. 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).
2. Winstock, A.R. and Mitcheson, L. New recreational drugs and the primary care approach to patients who use them. *BMJ* **344**, e288 (2012).
3. U. S. Department of Justice. GHB Analogs, *National Drug Intelligence Center* (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, 06/22/2023

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