

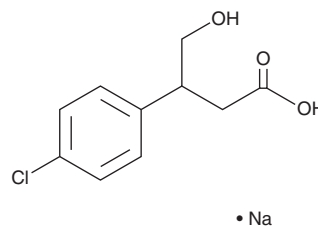
PRODUCT INFORMATION



3-(4-Chlorophenyl)-4-hydroxybutyric Acid (sodium salt)

Item No. 21483

CAS Registry No.: 1346603-21-1
Formal Name: 4-chloro-β-(hydroxymethyl)-benzenepropanoic acid, monosodium salt
Synonym: γ-Hydroxybaclofen
MF: C₁₀H₁₁ClO₃ • Na
FW: 237.6
Purity: ≥98%
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

3-(4-Chlorophenyl)-4-hydroxybutyric acid (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the 3-(4-chlorophenyl)-4-hydroxybutyric acid (sodium salt) in the solvent of choice, which should be purged with an inert gas. 3-(4-Chlorophenyl)-4-hydroxybutyric acid (sodium salt) is soluble in methanol. 3-(4-Chlorophenyl)-4-hydroxybutyric acid (sodium salt) is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

3-(4-Chlorophenyl)-4-hydroxybutyric acid is an inactive metabolite of the GABA_B receptor agonist baclofen (Item No. 18600).¹⁻³

References

1. Wuis, E.W., van Beijsterveldt, L.E., Dirks, R.J., *et al.* Rapid simultaneous determination of baclofen and its γ-hydroxy metabolite in urine by high-performance liquid chromatography with ultraviolet detection. *J. Chromatogr.* **420(1)**, 212-216 (1987).
2. Chéramy, A., Nieoullon, A., and Glowinski, J. Effects of the unilateral nigral application of baclofen on dopamine release in the two caudate nuclei of the cat. *Eur. J. Pharmacol.* **58(2)**, 133-140 (1979).
3. Ault, B. and Evans, R.H. The depressant action of baclofen on the isolated spinal cord of the neonatal rat. *Eur. J. Pharmacol.* **71(4)**, 357-364 (1981).

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, 10/10/2022

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