

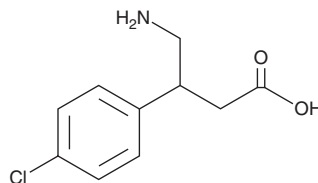
PRODUCT INFORMATION



(±)-Baclofen

Item No. 18600

CAS Registry No.: 1134-47-0
Formal Name: β-(aminomethyl)-4-chloro-benzenepropanoic acid
MF: C₁₀H₁₂ClNO₂
FW: 213.7
Purity: ≥98%
UV/Vis.: λ_{max}: 220, 266 nm
Supplied as: A crystalline 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

(±)-Baclofen is supplied as a crystalline solid. A stock solution may be made by dissolving the (±)-baclofen in water. The solubility of (±)-baclofen in water is approximately 4 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

(±)-Baclofen is a GABA_B receptor agonist (IC₅₀ = 180 nM).¹ It induces norepinephrine release from isolated rat atria (IC₅₀ = 4.5 μM). (±)-Baclofen (10 mg/kg) increases muscle rigidity in spastic mice.² It inhibits the norepinephrine-induced group II flexor reflex in anesthetized rats.³ (±)-Baclofen also reduces cocaine-induced hyperlocomotion in rats and binge-like ethanol intake in mice.^{4,5} Formulations containing baclofen have been used to treat muscle spasms caused by multiple sclerosis and spinal cord injury

References

1. Hill, D.R. and Bowery, N.G. ³H-baclofen and ³H-GABA bind to bicuculline-insensitive GABA_B sites in rat brain. *Nature* **290(5802)**, (1981).
2. Biscoe, T.J. and Fry, J.P. Some pharmacological studies on the spastic mouse. *Br. J. Pharmacol.* **75(1)**, 23-25 (1982).
3. Sakitani, K. The effects of centrally acting muscle relaxants on the intrathecal noradrenaline-induced facilitation of the flexor reflex mediated by group II afferent fibers in rats. *Jpn. J. Pharmacol.* **63(3)**, 369-376 (1993).
4. Lhuillier, L., Mombereau, C., Cryan, J.F., et al. GABA_B receptor-positive modulation decreases selective molecular and behavioral effects of cocaine. *Neuropsychopharmacology* **32(2)**, 388-398 (2007).
5. Moore, E.M. and Boehm, S.L., II Site-specific microinjection of baclofen into the anterior ventral tegmental area reduces binge-like ethanol intake in male C57BL/6J mice. *Behav. Neurosci.* **123(3)**, 555-563 (2009).

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, 11/28/2022

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