

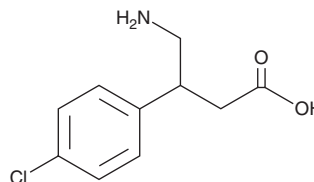
PRODUCT INFORMATION



(±)-Baclofen

Item No. 27326

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}: 258, 318, 332 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

(±)-Baclofen (Item No. 27326) is an analytical reference standard categorized as a gabapentinoid and a nootropic.^{1,2} (±)-Baclofen also has analgesic, skeletal muscle relaxant, and sedative properties.^{3,4} Formulations containing (±)-baclofen have been used in the treatment of spasticity in patients with multiple sclerosis or spinal cord injuries. This product is intended for use in analytical forensic applications. This product is also available as a general research tool (Item No. 18600).

References

1. Bowery, N.G., Hill, D.R., and Hudson, A.L. Characteristics of GABA_B receptor binding sites on rat whole brain synaptic membranes. *Br. J. Pharmacol.* **78(1)**, 191-206 (1983).
2. Tyurenkov, I.N., Borodkina, L.E., Bagmetova, V.V., et al. Comparison of nootropic and neuroprotective features of aryl-substituted analogs of gamma-aminobutyric acid. *Bull. Exp. Biol. Med.* **160(4)**, 465-469 (2016).
3. Chou, R., Peterson, K., and Helfand, M. Comparative efficacy and safety of skeletal muscle relaxants for spasticity and musculoskeletal conditions: A systematic review. *J. Pain Symptom Manage.* **28(2)**, 140-175 (2004).
4. de Beaurepaire, R., Sinclair, J.M.A., Heydtmann, M., et al. The use of baclofen as a treatment for alcohol use disorder: A clinical practice perspective. *Front. Psychiatry* **9**, 708 (2019).

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, 07/25/2023

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