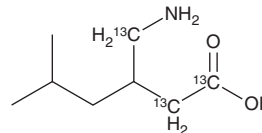


PRODUCT INFORMATION



(±)-Pregabalin-¹³C₃ Item No. 33879

CAS Registry No.: 1189980-48-0
Formal Name: 3-(aminomethyl-¹³C)-5-methylhexanoic-1,2-¹³C₂ acid
Synonyms: (±)-Isobutylgaba-¹³C₃, (R,S)-Isobutylgaba-¹³C₃
MF: C₅[¹³C]₃H₁₇NO₂
FW: 162.2
Purity: ≥98%
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.

Description

(±)-Pregabalin-¹³C₃ (Item No. 33879) is intended for use as an internal standard for the quantification for pregabalin (Item No. 13663) by GC- or LC-MS. Pregabalin is categorized as an anticonvulsant and analgesic.¹ It has potential for abuse and misuse.^{2,3} (±)-Pregabalin-¹³C₃ is regulated as a Schedule V compound in the United States. This product is intended for research and forensic applications.

References

1. Bryans, J.S. and Wustrow, D.J. 3-Substituted GABA analogs with central nervous system activity: A review. *Med. Res. Rev.* **19(2)**, 149-177 (1999).
2. Evoy, K.E., Morrison, M.D., and Saklad, S.R. Abuse and misuse of pregabalin and gabapentin. *Drugs* **77(4)**, 403-426 (2017).
3. Kapil, V., Green, J.L., Le Lait, M.C., et al. Misuse of the γ-aminobutyric acid analogues baclofen, gabapentin and pregabalin in the UK. *Br. J. Clin. Pharmacol.* **78(1)**, 190-191 (2014).

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, 04/11/2025

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