

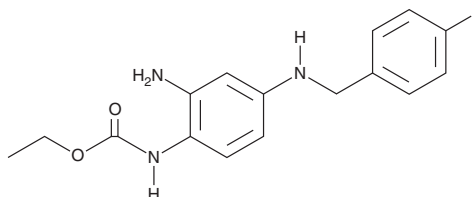
PRODUCT INFORMATION



Retigabine (exempt preparation)

Item No. 40314

CAS Registry No.: 150812-12-7
Formal Name: N-[2-amino-4-[[[4-(4-fluorophenyl)methyl]amino]phenyl]-carbamic acid, ethyl ester
Synonyms: D 23129, Ezogabine, GKE 841, GW 582892X, WAY-143841
MF: C₁₆H₁₈FN₃O₂
FW: 303.3
Purity: ≥98%
UV/Vis.: λ_{max}: 221, 254, 305 nm
Supplied as: A 1 mg/ml solution in acetonitrile
Storage: -20°C
Stability: ≥3 years



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

Description

Retigabine (exempt preparation) (Item No. 40314) is an analytical reference standard categorized as a positive allosteric modulator (PAM) of KCNQ2-5 (K_v7.2-7.5) potassium ion channels.¹ Formulations containing retigabine have been used in the treatment of partial epilepsies. Retigabine is regulated as a Schedule V compound in the United States. Retigabine (exempt preparation) (Item No. 40314) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

Reference

1. Gunthrope, M.J., Large, C.H., and Sankar, R. The mechanism of action of retigabine (ezogabine), a first-in-class K⁺ channel opener for the treatment of epilepsy. *Epilepsia* **53**(3), 412-424 (2012).

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, 01/15/2025

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