

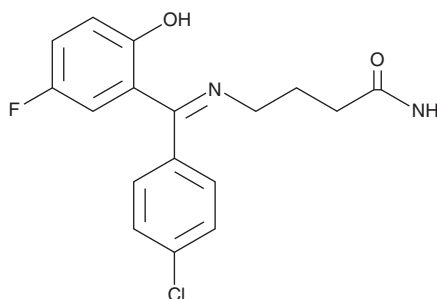
PRODUCT INFORMATION



Progabide

Item No. 45568

CAS Registry No.: 62666-20-0
Formal Name: 4-[[[(4-chlorophenyl)(5-fluoro-2-hydroxyphenyl)methylene]amino]butanamide
Synonym: SL 76002
MF: C₁₇H₁₆ClFN₂O₂
FW: 334.8
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

Progabide is supplied as a solid. A stock solution may be made by dissolving the Progabide in the solvent of choice, which should be purged with an inert gas. Progabide is sparingly soluble (1-10 mg/ml) in DMSO.

Description

Progabide is a GABA receptor agonist.¹ It selectively binds to the GABA_A and GABA_B receptors in postmortem rat cortical membranes (IC₅₀s = 35 and 100 μM, respectively) over α- and β-adrenergic, dopamine, serotonin, histamine H₁, and muscarinic receptors (IC₅₀s = >100 μM) in radioligand binding assays. Progabide (200 mg/kg) increases survival and reduces seizure progression to status epilepticus in rats.² It suppresses ethanol withdrawal-induced tremors and seizures in rats.³ Formulations containing progabide have been used in the treatment of epilepsy.

References

- Lloyd, K.G., Arbilla, S., Beaumont, K., *et al.* gamma-Aminobutyric acid (GABA) receptor stimulation. II. Specificity of progabide (SL 76002) and SL 75102 for the GABA receptor. *J Pharmacol Exp Ther.* **220**(3), 672-677 (1982).
- Mecarelli, O., de Feo, M.R., Rina, M.F., *et al.* Effects of progabide on bicuculline-induced epileptic seizures in developing rats. *Clin. Neuropharmacol.* **11**(5), 443-453 (1988).
- Fadda, F., Mosca, E., Meloni, R., *et al.* Suppression by progabide of ethanol withdrawal syndrome in rats. *Eur J. Pharmacol.* **109**(3), 321-325 (1985).

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, 05/13/2026

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