

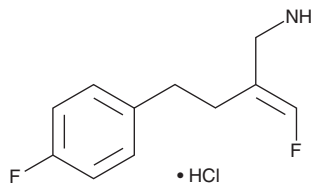
PRODUCT INFORMATION



Mofegiline (hydrochloride)

Item No. 35010

CAS Registry No.: 120635-25-8
Formal Name: 4-fluoro-βE-(fluoromethylene)-benzenebutanamine, monohydrochloride
Synonym: MDL 72974A
MF: C₁₁H₁₃F₂N • HCl
FW: 233.7
Purity: ≥95%
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

Mofegiline (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the mofegiline (hydrochloride) in water. The solubility of mofegiline (hydrochloride) in water is approximately 30 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Mofegiline is an inhibitor of monoamine oxidase B (MAO-B; IC₅₀ = 3.6 nM for the rat brain mitochondrial enzyme).¹ It is selective for MAO-B over MAO-A (IC₅₀ = 680 nM). Mofegiline also inhibits vascular adhesion protein-1 (VAP-1), also known as semicarbazide-sensitive amine oxidase (SSAO), an enzyme that has roles in leukocyte adhesion and transmigration, with an IC₅₀ value of 20 nM for the human enzyme.² It reduces MPTP-induced decreases in striatal levels of dopamine, dihydroxyphenylacetic acid (DOPAC), and homovanillic acid (HVA) in mice when administered at a dose of 1.25 mg/kg.¹ Mofegiline (5 mg/kg) also inhibits LPS-induced increases in bronchoalveolar lavage fluid (BALF) levels of TNF-α in transgenic mice overexpressing VAP-1.

References

1. Zreika, M., Fozard, J.R., Dudley, M.W., *et al.* MDL 72,974: A potent and selective enzyme-activated irreversible inhibitor of monoamine oxidase type B with potential for use in Parkinson's disease. *J. Neural Transm. Park. Dis. Dement. Sect.* **1(4)**, 243-254 (1989).
2. Foot, J.S., Deodhar, M., Turner, C.I., *et al.* The discovery and development of selective 3-fluoro-4-aryloxyallylamine inhibitors of the amine oxidase activity of semicarbazide-sensitive amine oxidase/vascular adhesion protein-1 (SSAO/VAP-1). *Bioorg. Med. Chem. Lett.* **22(12)**, 3935-3940 (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, 12/08/2022

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