

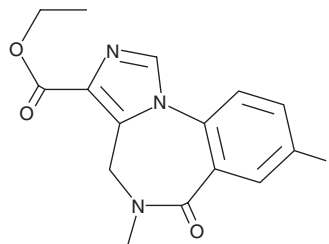
PRODUCT INFORMATION



Flumazenil

Item No. 14252

CAS Registry No.: 78755-81-4
Formal Name: 8-fluoro-5,6-dihydro-5-methyl-6-oxo-4H-imidazo[1,5-a][1,4]benzodiazepine-3-carboxylic acid, ethyl ester
Synonyms: Ro 1722, Ro 15-1788, Ro 41-8157
MF: C₁₅H₁₄FN₃O₃
FW: 303.3
Purity: ≥98%
UV/Vis.: λ_{max}: 244 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Flumazenil is supplied as a crystalline solid. A stock solution may be made by dissolving the flumazenil in the solvent of choice, which should be purged with an inert gas. Flumazenil is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of flumazenil in these solvents is approximately 1, 16, and 11 mg/ml, respectively.

Flumazenil is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, flumazenil should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Flumazenil has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Flumazenil is a GABA_A receptor antagonist (IC₅₀ = 2 nM in a radioligand binding assay using rat cortical synaptosomes).¹ Flumazenil also acts as a partial agonist of GABA_A receptors, decreasing the amplitude of electrically stimulated population spikes in rat hippocampal CA1 pyramidal neurons.² It increases the number of entries into the open arms of the elevated plus maze in high-anxiety BALB/c, but not C57BL/6, mice when administered at doses ranging from 0.1 to 1,000 μg/kg.³ Flumazenil (5 and 10 mg/kg) prevents a reduction in burying behavior induced by the GABA_A receptor positive allosteric modulator allopregnanolone (Item No. 16930) in ovariectomized rats when administered at doses of 5 and 10 mg/kg.⁴ Formulations containing flumazenil have been used to reverse sedation induced by benzodiazepines and in the treatment of benzodiazepine overdose or withdrawal.

References

1. Möhler, H., Burkard, W.P., Keller, H.H., *et al.* Benzodiazepine antagonist Ro 15-1788: Binding characteristics and interaction with drug-induced changes in dopamine turnover and cerebellar cGMP levels. *J. Neurochem.* **37(3)**, 714-722 (1981).
2. Buldalkova, S. and Weiss, M. Electrophysiological evidence for agonist properties of flumazenil, a benzodiazepine receptor antagonist, in rat hippocampus slices. *J. Neurol. Sci.* **149(2)**, 121-126 (1997).
3. Belzung, C., Le Guisquet, A.M., and Crestani, F. Flumazenil induces benzodiazepine partial agonist-like effects in BALB/c but not C57BL/6 mice. *Psychopharmacol. (Berl)*. **148(1)**, 24-32 (2000).
4. Fernández-Guasti, A. and Picazo, O. Flumazenil blocks the anxiolytic action of allopregnanolone. *Eur. J. Pharmacol.* **281(1)**, 113-115 (1995).

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, 01/17/2020

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