

Flumazenil

Benzodiazepine receptor ligand

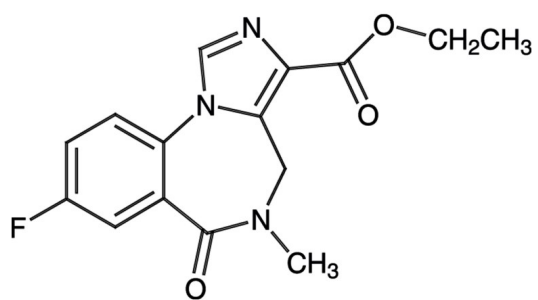
Benzodiazepine antagonist. Reverses all the effects of benzodiazepine agonists and inverse agonists.

Citations: 2

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Ro 15-1788
Appearance	White crystalline powder.
CAS	78755-81-4
Couple Target	Benzodiazepine receptor
Couple Type	Ligand
Formula	$C_{15}H_{14}FN_3O_3$
MW	303.3
Purity	≥98% (Assay)
Solubility	Soluble in ethanol (5 mg/ml warm) or DMSO (25 mg/ml warm)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com