

(+/-) Anatoxin

Nicotinic acetylcholine receptor ligand

(±)-Anatoxin is the synthetic version of (±)2-Acetyl-anatoxin, an alkaloidal toxin produced by the filamentous freshwater cyanophyte *Anabaena flos-aquae*. (±)-Anatoxin is a potent nicotinic agonist, stimulates dopamine release from rat striatal synaptosomes, and induces apoptosis in rat thymocytes and Vero cells following the generation of reactive oxygen species and the activation of caspases.

Citations: 7

[View Online »](#)

Ordering Information

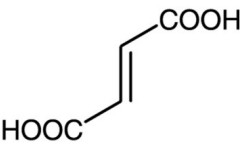
[Order Online »](#)

BML-C118-0001	1mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)

- Potent nicotinic agonist
- Toxin with apoptotic properties



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 1 month
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(±)-Anatoxin A fumarate, (±)-2-Acetyl-9-aza bicyclo[4,2,1]non-2-ene fumarate
Appearance	Light brown solid.
CAS	1219922-30-1
Couple Target	Acetylcholine receptor
Couple Type	Ligand
Formula	$C_{10}H_{15}NO \cdot C_4H_4O_4$
Identity	Identity determined by MS and ¹ H-NMR.
MW	281.3
Purity	≥96% (HPLC)
Solubility	Soluble in water (14mg/ml).
Source	Synthetic.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

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