

PRODUCT DATA SHEET



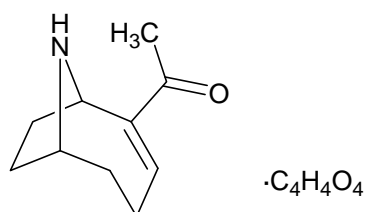
PRODUCT: (±)-Anatoxin A
2-Acetyl-9-aza-bicyclo[4.2.1]non-2-ene-fumarate

CAS NO. 64285-06-9

CATALOG NO.: C-118

LOT NO.: temp

STRUCTURE:



PHYSICAL APPEARANCE: White solid

MOLECULAR FORMULA: $C_{10}H_{15}NO \cdot C_4H_4O_4$

MOLECULAR WEIGHT: 281.3

PURITY: 98% (TLC)

SOLUBILITY: Soluble in water (14 mg/ml)

STORAGE: Store, as supplied, at $-20^{\circ}C$ for up to 1 year. Store solutions at $-20^{\circ}C$ for up to 3 months.

APPLICATION NOTES: An alkaloidal toxin produced by the filamentous freshwater cyanophyte *Anabaena flos-aquae*. It is a potent agonist at neuronal nicotinic acetylcholine receptors. At low concentrations (20-200 nM), it acts as an agonist. At higher concentrations (>800 nM), it blocks the AChR ion channel, desensitizes the receptor and blocks neuromuscular transmission.

NOTE: This vial may appear to be empty. This product is lyophilized. A clear, colorless film of anatoxin is deposited on the bottom of the vial.

- REFERENCES:
1. O.E.Edwards *et al. Can.J.Chem.* 1977 **55** 1367
 2. K.L.Swanson *et al. Mol.Pharmacol.* 1986 **29** 250
 3. P.Kofuji *et al. J.Pharmacol.Exp.Ther.* 1990 **252** 517
 4. P.Thomas *et al. J.Neurochem.* 1993 **60** 2308

The pharmacological and toxicological properties of this product have not been fully investigated. Exercise caution in use and handling. This product must not be used in humans.