

Acetylcholine receptor inhibitor; conjugated to tetramethylrhodamine

Instruction Manual

Catalog Number	PK-CA707-00012
Description	Red-fluorescent Tetramethylrhodamine- α -Bungarotoxin binds to the nicotinic acetylcholine receptor (AChR) with high affinity thereby inhibiting the binding of acetylcholine (ACh) to the receptor and blocking neuromuscular transmission. Thus, labeled alpha-bungarotoxin conjugates can be used to facilitate identification of nicotinic AChRs and image the distribution of the receptor at the neuromuscular junctions. See also our other dye- and biotin-labeled α -bungarotoxins, e.g. green-fluorescent Fluorescein- α -Bungarotoxin (PK-CA707-00011).
Quantity	500 μ g
Excitation / Emission Maxima	$\lambda_{ex}\backslash\lambda_{em} = 553/577$ nm
Molecular Structure	NA
Molecular Weight / Molecular Formula	~8500 Da; NA
Purity	NA
Appearance / Formulation / Solubility	Pink-red solid; soluble in water.
Storage & Stability	Store at -20°C. Protect from light, especially when in solution.
Applications	see Description
References	Anderson, et al., J. Cell Biol. 99, 1769(1984)
Caution	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.