

## Instruction Manual

Catalog Number	PK-CA707-00010-1
Description	Alpha-Bungarotoxin is isolated from the venom of the snake <i>Bungarus multicinctus</i> and is a potent inhibitor for the nicotinic acetylcholine receptors (motor endplate; $K_d = 1 \text{ nM}$ to $1 \text{ pM}$ ) of muscle cells at the neuromuscular junctions. The binding to the acetylcholine receptor sites is irreversible, producing neuromuscular blockade and skeletal muscle paralysis. It has become an important tool for the studies of neuromuscular junctions. See also our fluorescently labeled and biotin labeled bungarotoxins.
Quantity	1 mg
Excitation / Emission Maxima	NA
Molecular Structure	NA
Molecular Weight / Molecular Formula	7984 Da; $C_{338}H_{529}N_{97}O_{105}S_{11}$
Purity	Single band by FPLC.
Appearance / Formulation / Solubility	White solid; soluble in water.
Storage & Stability	Store desiccated at $-20^\circ\text{C}$ . Protect from light. Unstable at room temperature, especially when in solution.
Applications	see Description
References	Zhang, et al., <i>Neuron</i> , 12, 167(1994)
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.