

## Instruction Manual

Catalog Number	PK-CA707-00017
Description	Biotinylated $\alpha$ -bungarotoxin is useful in the affinity column isolation of the nicotinic AChR using an avidin or streptavidin agarose. Nicotinic AChRs labeled with biotin-XX- $\alpha$ -bungarotoxin can be localized using enzyme or fluorophore-labeled conjugates of avidin or streptavidin.
Quantity	500 $\mu$ g
Excitation / Emission Maxima	NA
Molecular Structure	NA
Molecular Weight / Molecular Formula	~8400 Da; NA
Purity	NA
Appearance / Formulation / Solubility	White lyophilized solid; soluble in water.
Storage & Stability	Store at -20°C. Protect from light, especially when in solution.
Applications	NA
References	1) J Cell Biol 125, 661-668 (1994) 2) J Biol Chem 268, 25108(1993) 3) Proc Natl Acad Sci U S A 77, 4823(1980)
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.