

Acetylcholine receptor inhibitor; conjugated to sulforhodamine 101
(also known as Texas Red*)

Instruction Manual

Catalog Number	PK-CA707-00015
Description	Sulforhodamine 101- α -bungarotoxin binds to the nicotinic acetylcholine receptor with high affinity. Since the dye sulforhodamine 101 has a longer absorption and emission wavelengths than tetramethylrhodamine- α -bungarotoxin (Cat.No. PK-CA707-00012), it gives a better spectral separation when a green fluorescent probe such as SynapseGreen C4 (FM1-43, Cat.No. PK-CA707-70020) is also used in the experiment. See also our other dye- and biotin-labeled α -bungarotoxins.
Quantity	500 μ g
Excitation / Emission Maxima	$\lambda_{ex}\backslash\lambda_{em} = 593/613$ nm
Molecular Structure	NA
Molecular Weight / Molecular Formula	~8600 Da; NA
Purity	NA
Appearance / Formulation / Solubility	Solid; soluble in water.
Storage & Stability	Store at -20°C. Protect from light, especially when in solution.
Applications	NA
References	1) J Cell Biol 150, 1385 (2000) 2) Neuron 23, 675(1999)
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.

* Texas Red is a trademark of Molecular Probes.

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