

## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** Biotin-XX- $\alpha$ -Bungarotoxin

**CATALOG #:** 00017

**MOLECULAR INFORMATION:** Mwt:  $\cong$  8400

**PROPERTIES:**

**Color & Form** White powder  
**Solubility** Soluble in water (dissolve in 0.5mL distilled water)  
**Absorption/Emission** N/A

**STORAGE AND HANDLING:** Store at -20 °C and protect from light, especially when in solution.

**APPLICATION:** Biotinylated  $\alpha$ -bungarotoxin is useful in the affinity column isolation of the nicotinic AChR using an avidin or streptavidin agarose.<sup>1</sup> Nicotinic AChRs labeled with biotin-XX- $\alpha$ -bungarotoxin can be localized using enzyme- or fluorophore-labeled conjugates of avidin or streptavidin.<sup>2,3</sup>

Ref.: 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).

**TOXICITY:** Not established.

<b>FIRST AID:</b>	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.
-------------------	---

<b>Disclaimer:</b> <i>Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.</i>
--