

PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Tetramethylrhodamine- α -Bungarotoxin

CATALOG #: 00012/00014

MOLECULAR INFORMATION: Mwt: \cong 8500

PROPERTIES:

Color & Form Pink red solid
Solubility Soluble in water
Absorption/Emission 553 nm \ 577 nm

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

APPLICATION: The labeled toxin binds to the nicotinic acetylcholine receptor with high affinity, and thus can be used to image the distribution of the receptor at the neuromuscular junctions.

Biotium also offers the green fluorescent fluorescein-alpha-bungarotoxin (**00011**).

Ref: Anderson, et al. *J. Cell Biol.* **99**, 1769(1984).

TOXICITY: Not established.

FIRST AID:	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.
-------------------	---

Disclaimer: <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>
--