

PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Sulforhodamine 101- α -Bungarotoxin

CATALOG #: 00015/00016

MOLECULAR INFORMATION: Mwt: \cong 8600

PROPERTIES:

Color & Form Bluish powder
Solubility Soluble in water
Absorption/Emission 593 nm / 613 nm

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

APPLICATION: 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 ($\lambda_{\text{ex}}/\lambda_{\text{em}} = 593/613$ nm) than tetramethylrhodamine- α -bungarotoxin (**00012**), it gives a better spectral separation when a green fluorescent probe such as SynptoGreen C4 (FM1-43, Cat# **70021**) is also used in the experiment.

Ref: 1) *J Cell Biol* **150**, 1385 (2000); 2) *Neuron* **23**, 675(1999)

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>
--