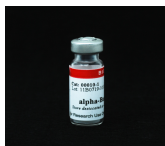


α -Bungarotoxin

A high affinity inhibitor of the nicotinic acetylcholine receptor in the neuromuscular junction. Blocks acetylcholine activity at the postsynaptic membrane.



Product Description

A potent neurotoxin that is an inhibitor for the motor endplate acetylcholine receptor ($K_d = 1 \text{ nM}$ to 1 pM). It binds at the neuromuscular junction with high affinity and is often used to study neuromuscular junctions by blocking cholinergic receptors. Alpha-bungarotoxin may also be used for detection of GABA A receptor subsets in cells (1), or for labeling recombinant proteins that express the alpha-bungarotoxin binding site (BBS) epitope tag (2).

- A polypeptide snake toxin
- Blocks and inhibits cholinergic receptors
- White solid soluble in water

Also see our [Fluorescent \$\alpha\$ -Bungarotoxins](#) and [Biotin \$\alpha\$ -Bungarotoxin](#) which can be used for labeling of nicotinic acetylcholine receptors at neuromuscular junctions in tissue sections. See our complete selection of α -Bungarotoxins below.

Product attributes

Toxin	Alpha-bungarotoxin
CAS number	11032-79-4
Molecular weight of antigen	7984
For live or fixed cells	For fixed cells, For live/intact cells
Cell permeability	Membrane impermeant
Storage Conditions	Store at -10 to -35 °C, Protect from light

α -Bungarotoxin, CF® Dye and Other Conjugates

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
Unconjugated	N/A	1 mg	00010-1	
Biotin-XX	N/A	0.5 mg	00017	
CF@405S	411/431 nm	100 ug	00002-100ug	CF@405S Features
0.5 mg	00002			
CF@488A	490/516 nm	100 ug	00005-100ug	CF@488A Features
0.5 mg	00005			
CF@543	543/563 nm	100 ug	00026-100ug	CF@543 Features
0.5 mg	00026			
CF@555	554/568 nm	100 ug	00018-100ug	CF@555 Features
0.5 mg	00018			
CF@568	562/584 nm	100 ug	00006-100ug	CF@568 Features
0.5 mg	00006			
CF@594	593/615 nm	100 ug	00007-100ug	CF@594 Features
0.5 mg	00007			
CF@633	629/650 nm	100 ug	00009-100ug	CF@633 Features
0.5 mg	00009			
CF@640R	642/663 nm	100 ug	00004-100ug	CF@640R Features
0.5 mg	00004			
CF@680R	680/701 nm	100 ug	00003-100ug	CF@680R Features
0.5 mg	00003			
Fluorescein (FITC)	498/517 nm	0.5 mg	00011	
10 x 50 ug	00013			
Tetramethylrhodamine (TRITC)	552/578 nm	0.5 mg	00012	
10 x 50 ug	00014			
Sulforhodamine-101 (Texas Red®)	595/613 nm	0.5 mg	00015	
10 x 50 ug	00016			

CF is a registered trademark of Biotium, Inc. Texas Red is a registered trademark of Thermo Fisher Scientific.

References

- PNAS, 103, 13, (2006), [DOI: 10.1073/pnas.0600847103](#)
- Meth. Enzymol., 521, (2013), [DOI: 10.1016/B978-0-12-391862-8.00006-5](#)
- Sci Adv, 6, 15, (2020), [DOI: 10.1126/sciadv.aax8382](#)
- Am J Physiol Cell Physiol, (2020) [DOI: 10.1152/ajpcell.00453.2019](#)
- Biochem Biophys Res Commun., 523, 214, (2020), [DOI: 10.1016/j.bbrc.2019.12.011](#)
- Front Cell Dev Biol., 8, 15, (2020), [DOI: 10.3389/fcell.2020.00015](#)
- Biophysics, 64, 772, (2019), [DOI: 10.1134/S0006350919050129](#)
- Cell Physiol Biochem, 53, 701, (2019), [DOI: 10.33594/000000166](#)
- Biochemistry (Mosc.), 4, 1085, (2019), [DOI: 10.1134/S0006297919090116](#)
- Biomaterials, 225, 119537, (2019), [DOI: 10.1016/j.biomaterials.2019.119537](#)
- ACS Chem. Biol., 13, 2568, (2018), [DOI: 10.1021/acschembio.8b00513](#)
- Cell Mol Neurobiol, 37, 1443, (2017), [DOI: 10.1007/s10571-017-0475-3](#)
- Neuroscience, 174, 234, (2011), [DOI: 10.1016/j.neuroscience.2010.11.016](#)
- Am J of Pathol, 177, 2509 (2010), [DOI: 10.2353/ajpath.2010.100243](#)
- Cell Tissue Biol, 4, 258, (2010), [DOI: 10.1134/S1990519X10030077](#)
- Neuroscience, 174, 234, (2009), [DOI: 10.1016/j.neuroscience.2010.11.016](#)
- J Cell Biol, 150, 1385, (2000), [DOI: 10.1083/jcb.150.6.1385](#)
- Neuron, 23, 675, (1999), [DOI: 10.1016/S0896-6273\(01\)80027-1](#)
- Neuron, 12, 167, (1994), [DOI: 10.1016/0896-6273\(94\)90161-9](#)
- J Cell Biol, 125, 661, (1994), [DOI: 10.1083/jcb.125.3.661](#)
- J Biol Chem, 268, 25108, (1993), [PMID: 8227074](#)
- Muscle Nerve, 5, 140, (1982), [DOI: 10.1002/mus.880050211](#)
- PNAS 77, 4823, (1980), [DOI: 10.1073/pnas.77.8.4823](#)
- Science, 196, 540, (1977), [DOI: 10.1126/science.850796](#)

This datasheet was generated on June 8, 2026 at 08:11:43 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/a-bungarotoxin-from-bungarus-multicinctus/>