

Molecular Probes™ α -Bungarotoxin and Conjugates User Guide

Catalog Numbers B56130, B13422, B35451, B13423, B35450, B1196, T1175, B1601

Pub. No. MAN0029177 Rev. A.0



WARNING! Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from [thermofisher.com/support](https://www.thermofisher.com/support).

Product Description

Extracted from *Bungarus multicinctus* venom, α -bungarotoxin has been used to stain acetylcholine receptors in skeletal muscle, rat myotubules, the electric organ from *Torpedo californica*, and transformed *Escherichia coli*.

Invitrogen™ offers unlabeled α -bungarotoxin and α -bungarotoxin labeled with biotin or a wide selection of fluorophores. (See Table 1)

Fluorescent α -bungarotoxin derivatives are useful for visualizing the distribution of acetylcholine receptors during neuromuscular junction development. The molecular weight of unlabeled α -bungarotoxin is about 8000 daltons. Approximately one fluorophore or one biotin is attached to each α -bungarotoxin, retaining optimal binding specificity. The labeled bungarotoxins are then chromatographically separated from unlabeled molecules to ensure maximum labeling of the product.

The Alexa Fluor™ 488 conjugate is the preferred green fluorescent derivative because of its brightness, pH sensitivity, and much greater photostability than fluorescein α -bungarotoxin. The far-red fluorescence of Alexa Fluor™ 647 and 680 conjugates permits three and four color experiments.

Contents and Storage

Table 1 α -Bungarotoxin and α -bungarotoxin conjugates

Material	Cat. No.	Amount ^[1]	Fluorescence excitation / emission	Storage
α -bungarotoxin, Alexa Fluor™ Plus 405 conjugate	B56130	100 μ g	411 nm / 431 nm	Store at $\leq 20^{\circ}\text{C}$. Desiccate. Avoid freeze-thaw cycles Protect from light
α -bungarotoxin, Alexa Fluor™ 488 conjugate	B13422	500 μ g	495 nm / 519 nm	
α -bungarotoxin, Alexa Fluor™ 555 conjugate	B35451	500 μ g	555 nm / 565 nm	
α -bungarotoxin, Alexa Fluor™ 594 conjugate	B13423	500 μ g	590 nm / 617 nm	
α -bungarotoxin, Alexa Fluor™ 647 conjugate	B35450	500 μ g	650 nm / 668 nm	
α -bungarotoxin, biotin-XX conjugate	B1196	500 μ g	N/A	
α -bungarotoxin, tetramethylrhodamine conjugate	T1175	500 μ g	554 nm / 577 nm	
α -bungarotoxin, unlabeled	B1601	1 mg (lyophilized)	N/A	

^[1] Unless otherwise noted, conjugates are lyophilized from phosphate-buffered saline, PBS, pH 7.4.

Required materials not supplied

Unless otherwise indicated, all materials are available through thermofisher.com. "MLS" indicates that the material is available from fisherscientific.com or another major laboratory supplier.

Catalog numbers that appear as links open the web pages for those products.

Item	Source
Phosphate-buffered saline, PBS, pH 7.4, 1 X 1000 mL	10010-031
Hank's balanced salt solution, HBSS, with no calcium chloride, magnesium chloride, magnesium sulfate, or phenol red, 1 x 1000 mL	14175-079
Image-iT™ Fixative Solution in PBS (4% formaldehyde, methanol-free), 1 X 20 mL	FB002
Image-iT™ Glyoxal 3% Fixative, 1 x 100 mL	I28700
Image-iT™ Paraformaldehyde 4% in 0.1 M Phosphate Buffer (Methanol free), 10 x 10 mL	I28800
Deionized water	MLS
Sodium	MLS
ProLong™ Glass Antifade Mountant, 5 x 2 mL	P36980b
EVOS™ M7000 Imaging System, 1	AMF7000

Prepare stock solution

- Prepare a 1.0 mg/mL α -bungarotoxin stock solution in PBS.
- The stock solution may be divided into aliquots and stored at -20°C for at least a month. For short term storage, add sodium azide to a final concentration of 2 mM, and store at 2-6°C.

Note: It is recommended that the protein conjugate solution is briefly centrifuged before use and that only the supernatant is added to the experiment. This step eliminates any protein aggregates that may have formed in the solutions, and reduces nonspecific background staining.

IMPORTANT! Protect from light and avoid repeated freezing and thawing.

Staining protocol

The following is an example protocol for staining 10 μ m-thick, fresh-frozen, cryosections of rat skeletal muscle with fluorescent α -bungarotoxin conjugates. This protocol may require optimization for other applications. The α -bungarotoxin conjugate staining can be performed concurrently with immunofluorescence staining.

1. Fix freshly frozen sections in 4% paraformaldehyde in PBS for 15 minutes at room temperature. Alternatively, sections can be fixed in ice-cold methanol for 5 minutes at -20°C. Rinse 3 times with PBS.
2. Permeabilize sections with PBS / 0.1% Triton-X™ 100 for 10 minutes at room temperature. Permeabilization is not required for methanol-fixed sections.
3. Prepare a 1 μ g/mL α -bungarotoxin staining solution in PBS. The conjugate can also be diluted in an immunofluorescence blocking buffer.
4. Rinse several times in PBS.
5. Mount in fluorescence antifade mounting medium and cover with a coverslip.
6. Rinse several times in PBS.
7. Mount in fluorescence antifade mounting medium and cover with a coverslip.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



Revision history: Pub. No. MAN0029177 A.0

Revision	Date	Description
A.0	1 March 2023	New document for Molecular Probes™ αConjugates.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2023 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.