

# $\alpha$ -Bungarotoxin and Conjugates

Rev. date 24 September 2012

Pub. no. MAN0006874

Detailed protocol is available online at [www.lifetechnologies.com/manuals](http://www.lifetechnologies.com/manuals).

**WARNING!** Before handling, read all applicable Safety Data Sheets (SDSs) at [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support).

- The molecular weight of unlabeled  $\alpha$ -bungarotoxin is approximately 8000.
- Approximately 1 fluorophore or 1 biotin is attached to each  $\alpha$ -bungarotoxin.
- Prepare a 1 mg/mL stock solution by reconstituting unlabeled  $\alpha$ -bungarotoxin in 1 mL of PBS or the  $\alpha$ -bungarotoxin conjugates in 0.5 mL of distilled water.

## Peak Excitation and Emission Wavelengths for $\alpha$ -Bungarotoxin Conjugates

Cat. No.	Conjugate	$\lambda_{\max}$ (nm)*	Em (nm)†
B1196	Biotin-XX	NA	NA
B13422	Alexa Fluor® 488	495	519
B13423	Alexa Fluor® 594	590	617
B35450	Alexa Fluor® 647	650	668
B35451	Alexa Fluor® 555	555	565
B35452	Alexa Fluor® 680	679	702
F1176	Fluorescein	494	518
T1175	Tetramethylrhodamine	554	577

\*Absorbance maximum for the dye. †Emission maximum for the dye.

### 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 found on Life Technologies' website at [www.lifetechnologies.com/termsandconditions](http://www.lifetechnologies.com/termsandconditions). If you have any questions, please contact Life Technologies at [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support).

© 2012 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners. LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

**For support visit** [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support)

[www.lifetechnologies.com](http://www.lifetechnologies.com)