

Streptavidin and Fluorescent Conjugates of Streptavidin

Table 1 Contents and storage

Material	Amount	Concentration	Storage	Stability
Unlabeled streptavidin*	5 mg	Lyophilized powder	• ≤-20°C • Desiccate	3 years
Fluorescent streptavidin conjugates	1 mg	Lyophilized powder	• ≤-20°C • Desiccate • Protect from light	6 months
	0.5 mL	2 mg/mL solution in PBS, pH 7.2, 5 mM sodium azide	• ≤6°C • Protect from light • Avoid freeze-thaw cycles	6 months
Streptavidin–R-PE and streptavidin–B-PE	1 mL	1 mg/mL solution in 0.1 M sodium phosphate, 0.1 M NaCl, 5 mM sodium azide, pH 7.5	2-8°CProtect from lightDo not freeze	6 months
Streptavidin-APC conjugates	0.5 mL (S868) or 250 µL (S32362)	1 mg/mL solution in 0.1 M sodium phosphate, 0.1 M NaCl, 5 mM sodium azide, pH 7.5	• 2-8°C • Protect from light • Do not freeze	6 months
Tandem conjugates of streptavidin	100 μL	1 mg/mL solution in 0.1 M sodium phosphate, 0.1 M NaCl, 2 mM EDTA, 1% glycerol, 5 mM sodium azide, pH 7.5	• 2–8°C • Protect from light • Do not freeze	6 months

^{*} The specific activity is >15 U/mg, where one unit is defined as the amount of protein required to bind 1 µg of biotin. Streptavidin is soluble to at least 10 mg/mL dissolved in phosphate-buffered saline (PBS) or other suitable buffer.

Introduction

Life Technologies carries an extensive line of unlabeled and labeled streptavidin products (Table 2, page 2), including conjugates of R-phycoerythrin (R-PE), B-phycoerythrin (B-PE), and allophyco-cyanin (APC) (Table 3, page 2). The high affinity of avidin, an egg-white protein, for biotin was first exploited in histochemical applications in the mid-1970s. ^{1,2} Avidin and its bacterial counterpart, streptavidin (from *Streptomyces avidinii*), have since become standard reagents for diverse detection schemes. ³ Streptavidin, a nonglycosylated 52,800-dalton protein, reportedly exhibits less nonspecific binding than avidin.

In addition to the products listed in this product information sheet, Life Technologies offers conjugates of avidin and NeutrAvidin® biotin-binding protein, fluorescent conjugates of many species-specific anti-IgG antibodies, and conjugates of protein A and protein G. For details, refer to our website at www.lifetechnologies.com or contact Technical Support.

For Research Use Only. Not for use in diagnostic procedures.

MAN0001787 | MP00888 Revision: 3.0

Table 2 Unlabeled streptavidin and fluorescent-dye conjugates of streptavidin

Label	Ex *	Em *	Catalog Number
Unlabeled streptavidin	NA	NA	S888
Fluorescent Dye Conjugates			
Alexa Fluor® 350	346	442	S11249
Marina Blue®	365	460	S11221
Cascade Yellow™	402	545	S11228
Alexa Fluor® 405	402	421	S32351
Pacific Blue™	410	455	S11222
Pacific Green™	411	500	S11200
Pacific Orange™	400	551	S32365
Alexa Fluor® 430	434	539	S11237
Fluorescein	494	518	S869
Alexa Fluor® 488	495	519	S11223, S32354
Alexa Fluor® 514	518	540	S32353
Alexa Fluor® 532	530	554	S11224
Alexa Fluor® 555	555	565	S21381, S32355
Tetramethylrhodamine	555	580	S870
Alexa Fluor® 546	556	573	S11225
Rhodamine B	570	590	S871
Rhodamine Red [™] -X	570	590	S6366
Alexa Fluor® 568	578	603	S11226
Alexa Fluor® 594	590	617	S11227, S32356
Texas Red®	595	615	S872
Texas Red®-X	595	615	S6370
Alexa Fluor® 633	632	647	S21375
Alexa Fluor® 635	633	647	S32364
Alexa Fluor® 647	650	668	S21374, S32357
Alexa Fluor® 660	663	690	S21377
Alexa Fluor® 680	679	702	S21378, S32358
Alexa Fluor® 700	702	723	S21383
Alexa Fluor® 750	749	775	S21384
Alexa Fluor® 790	784	814	S11378

^{*} Approximate fluorescence excitation (Ex) and emission (Em) maxima for conjugates, in nm. Complete spectra for most of these dyes are available at our website. † CMNB-caged fluorescein. NA = Not applicable.

Table 3 R-PE, B-PE, APC, and tandem conjugates of streptavidin

Label	Ex *	Em *	Catalog Number	
R-PE, B-PE, and Tandem-R-PE Conjugates				
B-Phycoerythrin (B-PE)	546, 565 †	575	S32350	
R-Phycoerythrin (R-PE)	496, 546, 565 †	578	S866, S21388 ‡	
Alexa Fluor® 610–R-PE	496, 546, 565 †	630	S20982	
Alexa Fluor® 647–R-PE	496, 546, 565 †	668	S20992	
APC and Tandem-APC Conjugates				
Allophycocyanin (APC)	650	660	S868, S32362 ‡	
Alexa Fluor® 700–APC	650	723	S21005	
Alexa Fluor® 750–APC	650	775	S21008	
**				

^{*} Approximate fluorescence excitation (Ex) and emission (Em) maxima for conjugates, in nm. † Multiple excitation peaks. ‡ Premium grade.

Unlabeled Streptavidin

Streptavidin has four binding sites, permitting a number of techniques in which unlabeled streptavidin can be used to bridge two biotinylated reagents. This bridging method, which is commonly used to link a biotinylated probe to a biotinylated enzyme in enzyme-linked immunohistochemical applications, often eliminates the background problems that can occur when using direct streptavidin–enzyme conjugates.

Fluorescent Streptavidin Conjugates

Some of the simplest fluorescent detection schemes involving streptavidin entail applying a biotinylated probe to the desired sample and then detecting the bound probe with a fluorescently labeled streptavidin. Fluorescent conjugates of streptavidin are commonly used to localize antigens in cells and tissues 4,5 and to detect biomolecules in immunoassays and DNA hybridization techniques.⁶⁻⁹ Peak excitation and emission wavelengths for these products are given in Table 2, page 2. The approximate degree of labeling of each conjugate is given in the certificate of analysis for each product lot.

R-PE, B-PE, APC and Tandem Conjugates of Streptavidin

Molecular Probes offers R-PE, B-PE, and APC conjugates of streptavidin, as well as tandem conjugates of R-PE and APC labeled with long-wavelength Alexa Fluor® dyes.

Our tandem conjugates (Table 3, page 2) comprise a donor phycobiliprotein, R-PE or APC, conjugated to longer-wavelength light-emitting fluorescence acceptors. By the process of fluorescence resonance energy transfer (FRET), an energy transfer cascade is established wherein most of the light absorbed by the donor R-PE or APC results in fluorescence of the acceptor dye. This process can be quite efficient, resulting in almost total transfer of energy to the acceptor dye.

Our premium grade R-PE and APC conjugates of streptavidin have been highly purified to ensure that the products are: 1) predominantly 1:1 conjugates of R-PE or APC and streptavidin; 2) free of all unconjugated streptavidin; and 3) mostly free of unconjugated R-PE or APC. The premium grade R-PE and APC conjugates are a fractionation of our S866 and S868 products, respectively, intended especially for applications that benefit from additional purity. The purity of each lot is confirmed by analytical HPLC.

The APC conjugates are prepared from chemically crosslinked APC to avoid dissociation of the molecule into subunits when highly diluted.¹⁰

Qdot® nanocrystal conjugates of streptavidin

For more information about our Qdot® nanocrystal streptavidin conjugates, visit www.lifetechnologies.com or contact Technical Support. The products are listed in Table 4, below.

Table 4 Qdot® conjugates of streptavidin

Label	Ex *	Em *	Catalog Number
Qdot® 525	<525	525	Q10143MP, Q10141MP
Qdot® 565	<565	565	Q10133MP, Q10131MP
Qdot® 585	<585	585	Q10113MP, Q10111MP
Qdot® 605	<605	605	Q10103MP, Q10101MP
Qdot® 625	<625	625	A10196
Qdot® 705	<705	705	Q10163MP, Q10161MP
Qdot® 800	<800	800	Q10173MP, Q10171MP

^{*} Qdot® nanocrystals are excitable (Ex, in nm) at any wavelength below their emission maxima (Em, in nm). For most practical applications, they should be exited at wavelengths below 450 nm. For additional information, refer to the product manual *Qdot*® *Streptavidin Conjugates*, available at www.lifetechnologies.com/manuals.

Preparing Unlabeled Streptavidin Conjugates

Reconstituted solutions are stable for approximately three months with the addition of sodium azide to a final concentration of 5 mM or thimerosal to 0.2 mM. For longer storage, divide solutions into aliquots and freeze at \leq -20°C. Avoid repeated freezing and thawing of solutions.

Preparing Labeled Streptavidin Conjugates

Solutions can be made by dissolving the powder in 0.5–1.0 mL of PBS or other suitable buffer. In the lyophilized powder form, the conjugates are stable for at least two years when stored as directed. Reconstituted solutions are stable for approximately six months with the addition of sodium azide to a final concentration of 5 mM or thimerosal to 0.2 mM. Fluorescent streptavidin conjugates supplied in solution are also stable for approximately six months. For longer storage, divide solutions into aliquots and freeze at <– 20° C. **Protect from light. Avoid repeated freezing and thawing of solutions.**

Labeling with Streptavidin Conjugates

Streptavidin conjugates are used as secondary detection reagents in histochemical applications, flow cytometry, ^{11,12} blot analysis, and immunoassays. These reagents can also be employed to localize biocytin, biocytin-X, biotin ethylenediamine and Alexa Fluor[®], Cascade Blue[®], or lucifer yellow biocytins—derivatives of biotin that are used as neuroanatomical tracers. ^{13,14} The following are commonly used methods for employing streptavidin as a secondary detection reagent.

- **Direct Procedure:** A biotinylated primary probe such as an antibody, single-stranded nucleic acid probe, or lectin is bound to tissues, cells, or other surfaces. Excess protein is removed by washing, and detection is mediated by reagents such as our fluorescent streptavidins or our enzyme-conjugated streptavidins plus substrate.
- Indirect Procedure: A biotinylated antibody or oligonucleotide is used to probe a tissue, cell, or other surface. This preparation is then treated with unlabeled streptavidin. Excess reagents are removed by washing, and detection is mediated by a biotinylated detection reagent such as fluorescein biotin (B1370), biotinylated R-PE (P811), biotinylated FluoSpheres® microspheres, or a biotinylated enzyme (P917) plus a substrate. Alternatively, an unlabeled primary antibody is bound to a cell followed by a biotinylated species-specific secondary antibody. After washing, the complex is detected by the direct or indirect procedures described above.

Centrifuge protein conjugate solutions briefly in a microcentrifuge before use and use only the supernatant for the experiment. This step will eliminate any protein aggregates that may have formed in solution, thereby reducing nonspecific background staining. Because staining protocols vary with the application, determine appropriate dilutions of conjugates empirically. For fluorescent dye conjugates of streptavidin, including R-PE, B-PE, APC, and tandem conjugates of streptavidin, a final concentration of 0.5–10 µg/mL is usually satisfactory for most histochemical applications.

Our Alexa Fluor® 680 and Alexa Fluor® 790 Streptavidin conjugates have been validated for use in western blot detection and are compatible with most standard fluorescent western scanners. These fluorescent markers can be used singly or in multicolor western applications. For membrane blot applications, Tween® 20 (0.1 to 0.2% v/v) and SDS (0.02% to 0.1% v/v) should be included in the detection buffer for the final incubation step to minimize non-specific background staining.

References

1. Proc Natl Acad Sci USA 71, 3537 (1974); 2. Biochim Biophys Acta 264, 165 (1972); 3. "Avidin-Biotin Technology," M. Wilchek and E. Bayer, Eds., Meth Enzymol 184 (complete volume); 4. J Cell Biol 111, 1183 (1990); 5. Physiol Plantarum 79, 231 (1990); 6. Cytometry 11, 126 (1990); 7. Proc Natl Acad Sci USA 87, 6223 (1990); 8. Science 249, 928 (1990); 9. Anal Biochem 171, 1 (1988); 10. Cytometry 8, 91 (1987); 11. J Microbial Methods 12, 1 (1990); 12. Biochemistry 16, 5150 (1977); 13. J Neurosci 10, 3421 (1990); 14. Brain Res 497, 361 (1989).

Product List Current prices may be obtained from our website or from our Customer Service Department.

Cat. no.	Product Name	Unit Size
S888	streptavidin	5 mg
S11249	streptavidin, Alexa Fluor® 350 conjugate	
S32351	streptavidin, Alexa Fluor® 405 conjugate	1 mg 1 mg
S11237	streptavidin, Alexa Fluor® 430 conjugate	
S11237 S11223	streptavidin, Alexa Fluor® 488 conjugate	1 mg
S32354	streptavidin, Alexa Fluor® 488 conjugate *2 mg/mL*	1 mg
		0.5 mL
S32352	streptavidin, Alexa Fluor® 500 conjugate	1 mg
S32353	streptavidin, Alexa Fluor® 514 conjugatestreptavidin, Alexa Fluor® 532 conjugate	1 mg
S11224		1 mg
S11225	streptavidin, Alexa Fluor® 546 conjugate	1 mg
S21381	streptavidin, Alexa Fluor® 555 conjugate	1 mg
S32355	streptavidin, Alexa Fluor® 555 conjugate *2 mg/mL*	0.5 mL
S11226	streptavidin, Alexa Fluor® 568 conjugate	1 mg
S11227	streptavidin, Alexa Fluor® 594 conjugate	1 mg
S32356	streptavidin, Alexa Fluor® 594 conjugate *2 mg/mL*	0.5 mL
S20982	streptavidin, Alexa Fluor® 610-R-phycoerythrin conjugate (Alexa Fluor® 610-R-phycoerythrin streptavidin) *1 mg/mL*	100 μL
S21375	streptavidin, Alexa Fluor® 633 conjugate	1 mg
S32364	streptavidin, Alexa Fluor® 635 conjugate	1 mg
S21374	streptavidin, Alexa Fluor® 647 conjugate	1 mg
S32357	streptavidin, Alexa Fluor® 647 conjugate *2 mg/mL*	0.5 mL
S20992	streptavidin, Alexa Fluor® 647-R-phycoerythrin conjugate (Alexa Fluor® 647-R-phycoerythrin streptavidin) *1 mg/mL*	100 μL
S21377	streptavidin, Alexa Fluor® 660 conjugate	1 mg
S21378	streptavidin, Alexa Fluor® 680 conjugate	1 mg
S32358	streptavidin, Alexa Fluor® 680 conjugate *2 mg/mL*	0.5 mL
S21383	streptavidin, Alexa Fluor® 700 conjugate	1 mg
S21005	streptavidin, Alexa Fluor® 700-allophycocyanin conjugate (Alexa Fluor® 700-allophycocyanin streptavidin) *1 mg/mL*	100 μL
S21384	streptavidin, Alexa Fluor® 750 conjugate	1 mg
S21008	streptavidin, Alexa Fluor® 750-allophycocyanin conjugate (Alexa Fluor® 750-allophycocyanin streptavidin) *1 mg/mL*	100 μL
S32363	streptavidin, Alexa Fluor® 750-R-phycoerythrin conjugate (Alexa Fluor® 750-R-phycoerythrin streptavidin) *1 mg/mL*	100 μL
S11378	streptavidin, Alexa Fluor® 790 conjugate	1 mg
S32362	streptavidin, allophycocyanin conjugate *premium grade* *1 mg/mL*	250 μL
S868	streptavidin, allophycocyanin, crosslinked, conjugate *1 mg/mL*	0.5 mL
S32350	streptavidin, B-phycoerythrin conjugate *1 mg/mL*	1 mL
S11228	streptavidin, Cascade Yellow™ conjugate	1 mg
S869 S11221	streptavidin, fluorescein conjugate	1 mg
	streptavidin, Marina Blue® conjugate	1 mg
S11222	streptavidin, Pacific Blue™ conjugate	1 mg
S11200	streptavidin, Pacific Green™ conjugate	500 μL
S32365	streptavidin, Pacific Orange™ conjugate	1 mg
S871	streptavidin, rhodamine B conjugate	1 mg
S6366	streptavidin, Rhodamine Red™-X conjugate	1 mg
S866	streptavidin, R-phycoerythrin conjugate *1 mg/mL*	1 mL
S21388	streptavidin, R-phycoerythrin conjugate *premium grade* *1 mg/mL*	1 mL
S870	streptavidin, tetramethylrhodamine conjugate	1 mg
S872	streptavidin, Texas Red® conjugate	1 mg
S6370	streptavidin, Texas Red®-X conjugate	1 mg

Purchaser Notification

Corporate Headquarters

5791 Van Allen Way Carlsbad, CA 92008

USA

Phone: +1 760 603 7200 Fax: +1 760 602 6500

Email: techsupport@lifetech.com

European Headquarters

Inchinnan Business Park 3 Fountain Drive Paisley PA4 9RF UK Phone: +44 141 814 6100

Toll-Free Phone: 0800 269 210 Toll-Free Tech: 0800 838 380 Fax: +44 141 814 6260 Tech Fax: +44 141 814 6117

Tech Fax: +44 141 814 6117 Email: euroinfo@invitrogen.com Email Tech: eurotech@invitrogen.com

Japanese Headquarters

LOOP-X Bldg. 6F 3-9-15, Kaigan Minato-ku, Tokyo 108-0022 Japan

Phone: +81 3 5730 6509 Fax: +81 3 5730 6519 Email: jpinfo@invitrogen.com

Additional international offices are listed at www.lifetechnologies.com

These high-quality reagents and materials must be used by, or directly under the supervision of, a technically qualified individual experienced in handling potentially hazardous chemicals. Read the Safety Data Sheet provided for each product; other regulatory considerations may apply.

Obtaining Support

For the latest services and support information for all locations, go to www.lifetechnologies.com.

At the website, you can:

- Access worldwide telephone and fax numbers to contact Technical Support and Sales facilities
- Search through frequently asked questions (FAQs)
- Submit a question directly to Technical Support (techsupport@lifetech.com)
- Search for user documents, SDSs, vector maps and sequences, application notes, formulations, handbooks, certificates of
 analysis, citations, and other product support documents
- · Obtain information about customer training
- · Download software updates and patches

SDS

Safety Data Sheets (SDSs) are available at www.lifetechnologies.com/sds.

Certificate of Analysis

The Certificate of Analysis provides detailed quality control and product qualification information for each product. Certificates of Analysis are available on our website. Go to www.lifetechnologies.com/support and search for the Certificate of Analysis by product lot number, which is printed on the product packaging (tube, pouch, or box).

Disclaimer

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.

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. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

Limited Use Label License: Research Use Only

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial services of any kind, including, without limitation, reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact **outlicensing@lifetech.com** or Out Licensing, Life Technologies Corporation, 5791 Van Allen Way, Carlsbad, California 92008.

The trademarks mentioned herein are the property of Life Technologies Corporation and/or its affiliate(s) or their respective owners.

Cy is a registered trademark of GE Healthcare Limited UK.
NeutrAvidin is a registered trademark of Pierce Biotechnology, Inc.
Tween is a registered trademark of Uniqema Americas LLC.

©2013 Life Technologies Corporation. All rights reserved.

