

Human Neural Stem Cell Immunocytochemistry Kit

Catalog no. A24354

Table 1 Contents and storage

Kit component	Part no.	Concentration	Amount	Storage	Usage notes
Primary antibodies					
anti-NESTIN (host: mouse)	A24345	- - 50X	20 μL	-20°C to 4°C	Dilute with Blocking Solution
anti-PAX6 (host: rabbit)	A24340				
anti-SOX1 (host: goat)	A24347				
anti-SOX2 (host: rabbit)	A24339				
Secondary antibodies					
Alexa Fluor® 488 donkey anti-mouse; for use with anti-NESTIN	A24350	250X	20 μL	–20°C to 4°C; avoid freeze-thaw cycles	Ex/Em ¹ 495/519 nm (green); spin before use ²
Alexa Fluor® 488 donkey anti-goat; for use with anti-SOX1	A24349				Ex/Em ¹ 495/519 nm (green); spin before use ²
Alexa Fluor® 555 donkey anti-rabbit; for use with anti-PAX6 or anti-SOX2	A24342		40 μL		Ex/Em ¹ 555/565 nm (orange); spin before use ²
Alexa Fluor® 594 donkey anti-rabbit; for use with anti-PAX6 or anti-SOX2	A24343				Ex/Em ¹ 590/617 nm (red); spin before use ²
Additional reagents					
NucBlue [®] Fixed Cell Stain (DAPI nuclear DNA stain)	R37606	NA	1 vial	–20°C to ambient temperature	Ex/Em ¹ 358/461 nm (blue); apply 2 drops/mL
Fixative Solution	A24344	1X	10 mL		4% formaldehyde in DPBS
Permeabilization Solution	A24352				0.5% Triton® X-100 in DPBS
Blocking Solution	A24353		20 mL		3% BSA in DPBS
Wash Buffer	A24348	10X			Dilute to 1X with water

Handling and shelf life: Use aseptic technique when handling all reagents. Allow frozen reagents to thaw completely before using them. Once thawed, the kit should not be refrozen (aliquots not recommended). Store at 4°C for up to 6 months.

Description

This kit contains a complete set of primary and secondary antibodies along with ready-to-use buffers to enable facile immunocytochemistry characterization of human neural stem cells (NSC). The primary antibodies included in this kit target well-established NSC markers (Nestin, PAX6, SOX1, and SOX2) and were carefully selected to help ensure excellent performance in immunocytochemistry applications.

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

MAN009671 Revision A.0

¹ Approximate excitation/emission wavelength maxima.

 $^{^2}$ Centrifuge Secondary Antibody solutions (e.g., 2 minutes at 10,000 \times g) and add only the supernatant to the Blocking Solution. This step eliminates any protein aggregates that may have formed during storage, thereby reducing non-specific background staining.

See Table 2, page 3, for recommended volumes to use based on the culture format of the cells to be stained. See Table 3, page 3, for multiplex staining options.

Caution: Use gentle liquid handling and pipeting techniques when adding or removing liquids to minimize the possibility of dislodging cells and losing them during the handling steps.



1. Remove media from the cells.





2. Add Fixative Solution and incubate for 15 minutes at room temperature.



3. Remove Fixative Solution.

Optional stopping point: After removing Fixative, add Wash Buffer (diluted to 1X with water), wrap the sample to prevent it from drying out, and store at 4°C for up to 2 weeks.





Add Permeabilization Solution and incubate 15 minutes at room temperature.



5. Remove Permeabilization Solution.





6. Add Blocking Solution and incubate 1 hour at room temperature.



7. Remove Blocking Solution.





Add desired Primary Antibody (diluted to 1X in Blocking Solution; see Table 3 for options) and incubate for 3 hours at room temperature (or overnight at 4°C).



9. Wash 3 times with Wash Buffer (diluted to 1X with water).





10. Add the appropriate Secondary Antibody (diluted to 1X in Blocking Solution; see Table 3 for guidance) and incubate for 1 hour at room temperature.



11. Wash 3 times with Wash Buffer (diluted to 1X with water).

Optional: Add 2 drops/mL of NucBlue® Fixed Cell Stain (DAPI) into the last wash step and incubate for 5 minutes.



12. Image the cells immediately. Cells can also be stored at 4°C in the dark for up to 1 week.

Table 2 Recommended final volumes to use during the protocol.

Culture format	Volume	No. of tests ¹
96-well plate	50 μL/well	20
48-well plate	100 μL/well	10
24-well plate	200 μL/well	5
12-well plate	400 μL/well	2
6-well plate	1000 μL/well	1
35-mm dish	1000 μL/dish	1
4-well chamber slide	400 μL/well	2
8-well chamber slide	200 μL/well	5

¹ When using the suggested staining volume, this kit contains sufficient reagents for the indicated number of tests per primary antibody.

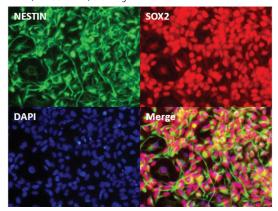
Table 3 Dual antibody staining options. Note that the NucBlue® Fixed Cell Stain (a DAPI nuclear DNA stain) provided in this kit is also compatible with these antibody combinations. See Figure 1, below, for example pictures.

Color options:	Green ¹ (e.g., FITC filter)	Orange ¹ (e.g., Cy [®] 3 / TRITC filter) or Red ¹ (e.g., Texas Red [®] filter)			
Antibody combination # 1: NESTIN + SOX2					
Primary antibody	anti- NESTIN (host: mouse)	anti-SOX2 (host: rabbit)			
Secondary antibody	Alexa Fluor® 488 donkey anti-mouse	Alexa Fluor [®] 555 donkey anti-rabbit or Alexa Fluor [®] 594 donkey anti-rabbit			
Antibody combination # 2: SOX1 + PAX6					
Primary antibody	anti-SOX1 (host: goat)	anti- PAX6 (host: rabbit)			
Secondary antibody	Alexa Fluor® 488 donkey anti-goat	Alexa Fluor [®] 555 donkey anti-rabbit or Alexa Fluor [®] 594 donkey anti-rabbit			
¹ See Table 1, page 1, for	approximate excitation/emission wavelength	maxima.			

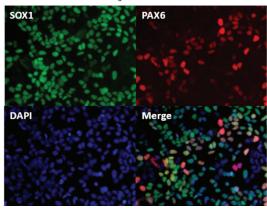
Figure 1 Neural stem cells derived from an iPSC line using Gibco® PSC Neural Induction Medium (Cat. no. A1647801) were stained for NSC markers Nestin and SOX2 (antibody combination #1) or SOX1 and PAX6 (antibody combination #2) and nuclear DNA (DAPI) using the Human Neural Stem Cell Immunocytochemistry Kit (Cat. no. A24354).

Note that not all NSCs will stain positive for PAX6, this is normal and to be expected (i.e., only NSCs that possess forebrain neuron fate commitment will stain positive). In our experience, approximately 15–50% of NSCs generated using Gibco® PSC Neural Induction Medium stain positive for PAX6.

Antibody combination # 1: NESTIN + SOX2 with additional DAPI (nuclear DNA) staining.



Antibody combination # 2: SOX1 + PAX6 with additional DAPI (nuclear DNA) staining



Cat. no.	Product Name	Unit Size
A24354	Human Neural Stem Cell Immunocytochemistry Kit	1 kit
Related Pro	ducts	
A1647801	Gibco® PSC Neural Induction Medium	1 kit
N7800100	Gibco® Human Neural Stem Cells (H9-Derived)	1 mL
R37602	Image-iT® Fixation/Permeabilization Kit	1 kit
A13998	OCT4 Rabbit Monoclonal Antibody (clone C30A3)*	100 μL
	*Pluripotent marker OCT4 is a good negative control marker for NSCs	

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

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
Temail: 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).

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.

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.

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.

The trademarks mentioned herein are the property of Life Technologies Corporation and/or its affiliate(s) or their respective owners. Triton is a registered trademark of Union Carbide Corporation.

Cy3 is a registered trademarks of GE Healthcare UK, Ltd.

©2013 Life Technologies Corporation. All rights reserved.

