

True-Nuclear™ Human Treg Flow™ Kit (FOXP3 Alexa Fluor® 488/CD4 PE-Cy5/CD25 PE)

Catalog# / Size	320027 / 25 tests
Clone	150D
Regulatory Status	RUO
Other Names	Forkhead box protein P3, Scurfin, JM2, IPEX, Zinc finger protein JM2

Description T regulatory (Treg) cells are a subset of T lymphocytes which is characterized by CD4+/CD25+/FOXP3+. These naturally occurring Treg cells originate in the thymus, and comprise 2-10% of peripheral CD4+ T cells. It has been shown that Treg cells are able to inhibit T cells proliferation and cytokine production and play critical roles in preventing autoimmunity as well as in controlling tumor immunity and transplantation tolerance. Impaired Treg function or Treg cell deficiency will develop variety of autoimmune diseases, while higher frequency of Treg cells will cause hypo-immune response to pathogens.

BioLegend's True-Nuclear™ Human Treg Flow™ Kit is designed and formulated specifically for immunofluorescence staining and flow cytometric analysis of human Treg cells in a mixed lymphocyte population. This kit is composed of fluorochrome conjugated anti-human CD4, CD25, FOXP3 antibodies, and the critical buffers. It is easy to use for identification of Treg cells.

Product Details

Verified Reactivity	Human
Antibody Type	Monoclonal
Concentration	Lot-specific (to obtain lot-specific concentration, please enter the lot number in our Concentration and Expiration Lookup or Certificate of Analysis online tools.)
Storage & Handling	This kit is guaranteed for six months. Upon receipt, store between 2°C and 8°C, and protected from prolonged exposure to light. Do not freeze.
Application	ICFC - Quality tested
Application Notes	<p>Materials Provided:</p> <ol style="list-style-type: none">1. Alexa Fluor® 488 anti-mouse/rat/human FOXP3 - 25 tests2. Alexa Fluor® 488 Mouse IgG1, κ isotype control - 25 tests3. True-Nuclear™ Transcription Factor Buffer Set - 120 tests (Cat. No. 424401)4. anti-human CD4 PE-Cy5/CD25 PE Cocktail - 50 tests <p>Materials Not Included:</p> <ol style="list-style-type: none">1. Cell Staining Buffer (Cat. No. 420201)2. Single color compensation controls <p>Immunofluorescence Staining Procedures:</p> <ol style="list-style-type: none">1. Perform cell surface staining as described in BioLegend's Cell Surface Immunofluorescence Staining Protocol. Add 20 µL of the anti-human CD4 PE-Cy5/CD25 PE cocktail to each tube and incubate in the dark for 20 minutes.2. Add 2 mL of the cell staining buffer, centrifuge tubes at 400 x g at room temperature for five minutes, and discard the supernatant.3. Repeat Step 2, for a total of two washes.4. Add 1 mL of the Transcription Factor 1X Fix solution to each tube, vortex, and incubate at room temperature in the dark for 45-60 minutes.5. Without washing, add 2 mL of the Transcription Factor 1X Perm Buffer to each tube.6. Centrifuge tubes at 400 x g at room temperature for five minutes, and discard the supernatant.7. Add 2 mL of the Transcription Factor 1X Perm Buffer to each tube.8. Centrifuge tubes at 400 x g at room temperature for five minutes, and discard the supernatant.9. Resuspend the cell pellet in 100 µL of the Transcription Factor 1X Perm Buffer.10. Add 5 µL of Alexa Fluor® 488 anti-mouse/rat/human FOXP3 antibody or 5 µL of Alexa Fluor® 488 mouse IgG1, κ isotype control into the appropriate tubes. Incubate in the dark at room temperature for at least 30 minutes.11. Add 2 mL of the Transcription Factor 1X Perm Buffer to each tube.

12. Centrifuge tubes at 400 x g at room temperature for five minutes, and discard the supernatant.
13. Add 2 mL of the cell staining buffer.
14. Centrifuge tubes at 400 x g at room temperature for five minutes, and discard the supernatant.
15. Resuspend in 0.5 mL cell staining buffer and then acquire tubes on a flow cytometer.

Caution: The True-Nuclear™ Transcription Factor Buffer Set contains paraformaldehyde, which is toxic and mutagenic. Please handle with caution. Wear gloves, lab coats, and necessary protection to avoid direct contact.

NOTE: For flow cytometric staining with this clone, True-Nuclear™ Transcription Factor Buffer Set (Cat. No. [424401](#)) offers improved staining and is highly recommended over the Foxp3/Perm Buffer Set (Cat. No. [421403](#)).

Application References

(PubMed link indicates BioLegend citation)

1. Roncador G, et al. 2005 *Eur. J. Immunol.* 35:1681.
2. Mayack. S, et al. 2006. *J. Immunol.* 176:2059. [J. Immunol](#)
3. Yang ZZ, et al. 2006. *Blood* 107:3639.
4. Gavin MA, et al. 2006. *P. Natl. Acad. Sci. USA* 103:6659.
5. Groh V, et al. 2006. *Nature Immunology* 7:755.
6. Lewkowicz P, et al. 2006 *J. Immunol.* 177:7155.
7. Luke PPW, et al. 2006. *Amer. J. Transplant.* 6(9):2023.
8. Bamias G, et al. 2007. *J. Immunol.* 178:1809.
9. Valencia X, et al. 2007. *J. Immunol.* 178:2579. [PubMed](#)
10. Davidson TS, et al. 2007. *J. Immunol.* 178:4022.
11. MacDonald K PA, et al. 2007. *Blood* doi:10.1182/blood-2007-01-067249.
12. Jaffar Z, et al. 2007. *J. Immunol.* 179:6193.
13. Müller M, et al. 2007. *J. Immunol.* 179:2774.
14. Jordan JM, et al. 2008. *Infect Human.* 76:3717. [PubMed](#)
15. Golovina TN, et al. 2008. *J. Immunol.* 181:2855. [PubMed](#)
16. Fallarino F, et al. 2009. *J. Exp Med.* 206:2511. [PubMed](#)
17. Banham Alison, et al. 2009. *Vet Immunol and Immunop* 127.3-4:376-381
18. Klunker S, et al. 2009. *J. Exp Med.* [PubMed](#)
19. Haque A, et al. 2010. *J. Immunol.* 184:2583. [PubMed](#)
20. Liu Y, et al. 2012. *Food Chem Toxicol.* 50:1920. [PubMed](#)

"

Antigen Details

Distribution	T regulatory cells
Cell Type	Tregs
Biology Area	Cell Biology, Immunology, Transcription Factors
Molecular Family	CD Molecules, Nuclear Markers
Antigen References	<ol style="list-style-type: none"> 1. Hori S, et al. 2003. <i>Science</i> 299:1057. 2. Fontenot JD, et al. 2003. <i>Nature Immunol.</i> 4:330. 3. Ferguson PJ, et al. 2000. <i>Am. J. Med. Genet.</i> 90:390. 4. Bennett CL, et al. 2001. <i>Nature Genet.</i> 27:20. 5. Allan SE, et al. 2005. <i>J. Clin. Invest.</i> 115:3276.

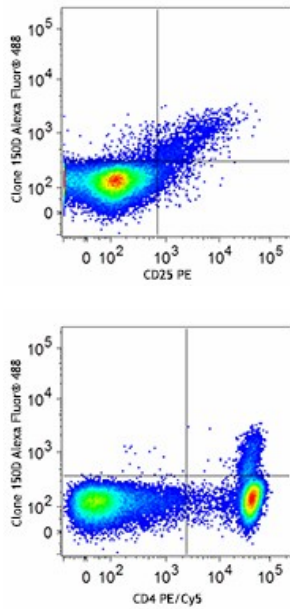
Gene ID	20371
	317382
	50943

Other Formats

Purified anti-mouse/rat/human FOXP3, PE anti-mouse/rat/human FOXP3, Alexa Fluor® 488 anti-mouse/rat/human FOXP3, Alexa Fluor® 647 anti-mouse/rat/human FOXP3, True-Nuclear™ Human Treg Flow™ Kit (FOXP3 Alexa Fluor® 488/CD4 PE-Cy5/CD25 PE), True-Nuclear™ Mouse Treg Flow™ Kit (FOXP3 Alexa Fluor® 488/CD4 APC/CD25 PE)

Product Data

Human peripheral blood lymphocytes were stained with True-Nuclear™ Human Treg Flow™ Kit (FOXP3 Alexa Fluor® 488/CD4 PE-Cy5/CD25 PE).



For research use only. Not for diagnostic use. Not for resale. BioLegend will not be held responsible for patent infringement or other violations that may occur with the use of our products.

*These products may be covered by one or more Limited Use Label Licenses (see the BioLegend Catalog or our website, www.biolegend.com/ordering#license). BioLegend products may not be transferred to third parties, resold, modified for resale, or used to manufacture commercial products, reverse engineer functionally similar materials, or to provide a service to third parties without written approval of BioLegend. By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.

BioLegend Inc., 8999 BioLegend Way, San Diego, CA 92121 www.biolegend.com
Toll-Free Phone: 1-877-Bio-Legend (246-5343) Phone: (858) 768-5800 Fax: (877) 455-9587