

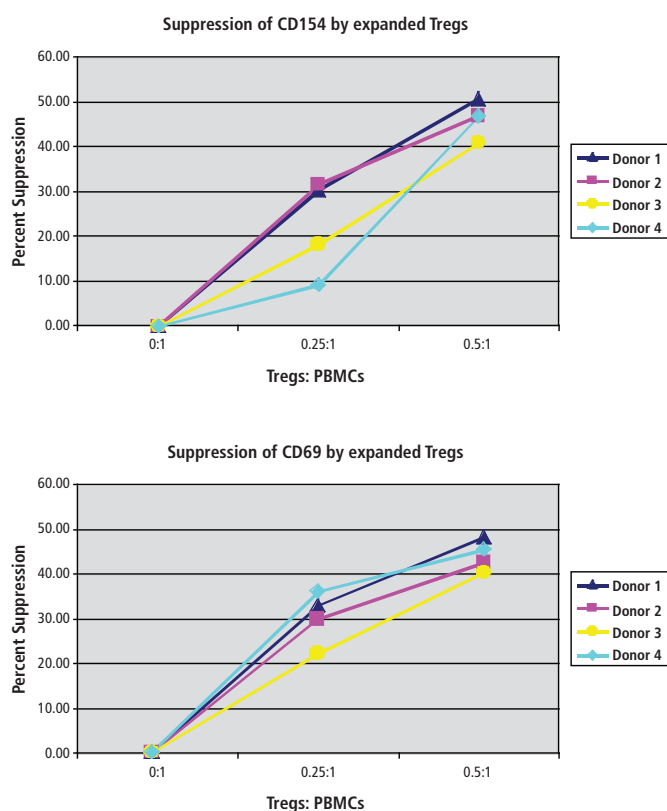
# BD FastImmune Human Regulatory T Cell Function Kit

## Features

Assess the function of expanded regulatory T cells (Tregs) in just one day

Detects the Treg-mediated suppression of the activation markers CD69 and CD154 on responding T cells using flow cytometry

Provided as a convenient cocktail of CD3, CD4, and CD25 with CD154 and CD69 as drop-in reagents



**Figure 1.** Treg-mediated suppression of CD154 (top) and CD69 (bottom) (n=4).

CD4<sup>+</sup>CD25<sup>high</sup>CD127<sup>low</sup>CD45RA<sup>+</sup> sorted Tregs were expanded in culture for 14 days. At a 0.5:1 Treg:PBMC ratio, Tregs suppressed the frequency of CD4<sup>+</sup>CD154<sup>+</sup> and CD4<sup>+</sup>CD69<sup>+</sup> lymphocytes by approximately 46% and 44% (on average) respectively, when PBMCs were stimulated with anti-CD3/CD28 beads for 7 hours.

As published data on the immunosuppressive potential of Regulatory T Cells (Tregs) has accumulated, enthusiasm for their potential application has intensified. BD Biosciences has a large portfolio of products relevant to Treg research, including this BD FastImmune™ Human Regulatory T Cell Function Kit, and the portfolio continues to expand to support new discoveries.

Tregs are a rare cell population representing 5–10% of all CD4<sup>+</sup> cells, and prevent aberrant immune responses. Low numbers of Tregs, or functionally defective Tregs, can lead to autoimmune diseases such as type I diabetes and graft-vs-host disease (GvHD). Tregs suppress effector T cells by inhibiting activation, proliferation, differentiation, and effector function.<sup>1,2</sup>

The Human Regulatory T Cell Function Kit measures the expression of activation markers CD69 and CD154 on effector T cells. A decrease in the levels of activation markers in the presence of Tregs may indicate the suppressive function of Tregs.

## How the assay works

Varying concentrations of expanded Tregs are co-cultured with responding effector cells in the presence of T-cell specific stimuli such as CD3/CD28 beads. To capture transiently expressed CD154 on the cell surface, APC-labeled CD154 reagent is present during the entire stimulation time. After seven hours of activation, the frequency of CD154-positive and/or CD69-positive effector T cells is compared in the presence and absence of Tregs.

The BD FastImmune Human Regulatory T Cell Function Kit contains a 3-color cocktail of CD3 PerCP-Cy™5.5 (clone SK7), CD4 FITC (clone SK3), CD25 PE (clone 2A3), and single color CD154 APC (clone 89-76) and CD69 PE-Cy™7 (clone L78) as drop-in reagents. BD FastImmune EDTA is also included. To accommodate various flow cytometer configurations, this assay can be stained as a 4-color (by excluding CD69 PE-Cy7) or a 5-color experiment.

## Tools for the sorting of Tregs

In addition to the Treg functional assay, the BD Human Regulatory T Cell Sorting Kit is also available. The kit contains a 3-color reagent cocktail that can be used to identify and sort CD4<sup>+</sup>CD25<sup>int/high</sup>CD127<sup>low</sup> expandable natural Treg populations. The reagent cocktail includes CD4 PerCP-Cy5.5 (clone L200), CD25 PE (clone 2A3), and CD127 Alexa Fluor® 647 (clone 40131). The one-step pre-mixed cocktail simplifies Treg identification and significantly enhances enrichment of Treg populations by 2 to 4 times compared to gating on CD25<sup>high</sup> alone.

This kit also includes CD45RA FITC (clone HI100) as an optimized drop-in. Publications by Hoffmann<sup>3</sup> and Miyara<sup>4</sup> have reported that isolation of the CD45RA<sup>+</sup> subset of Tregs may provide an increased percentage of FoxP3<sup>+</sup> Tregs with greater suppressive capacity post enrichment compared to CD45<sup>-</sup> Tregs.

Visit [bdbiosciences.com/tcell](http://bdbiosciences.com/tcell) for more information.

# BD FastImmune Human Regulatory T Cell Function Kit

## References

1. Sakaguchi S, Wing K, Onishi Y, Prieto-Martin P, Yamaguchi T. Regulatory T cells: how do they suppress immune responses? *Int Immunol*. 2009;21:1105-1111.
2. Askenasy N, Kaminitz A, Yarkoni S. Mechanisms of T regulatory cell function. *Autoimmun Rev*. 2008;7:370-375.
3. Hoffmann P, Eder R, Boeld TJ, et al. Only the CD45RA+ subpopulation of CD4+CD25high T cells gives rise to homogeneous regulatory T-cell lines upon in vitro expansion. *Blood*. 2006;108:4260-4267.
4. Miyara M, Yoshioka Y, Kitoh A, et al. Functional delineation and differentiation dynamics of human CD4+ T cells expressing the FoxP3 transcription factor. *Immunity*. 2009;30:899-911.

## Ordering Information

Description	Contents	Size	Cat. No.
Human Regulatory T Cell Function Kit	Human Regulatory T Cell Function Cocktail (CD3 PerCP-Cy5.5, CD4 FITC, CD25 PE), CD154 APC, CD69 PE-Cy7, and EDTA Solution	50 Tests	648956

## Related Products

Description	Contents	Size	Cat. No.
Human Regulatory T Cell Sorting Kit	Human Regulatory T Cell Sorting Cocktail (CD4 PerCP-Cy5.5, CD25 PE, CD127 Alexa Fluor® 647), CD45RA FITC, and compensation controls	10 Tests 1 Test = 1 x 10 <sup>6</sup> cells	560753



Alexa Fluor® is a registered trademark of Molecular Probes, Inc.

Cy™ is a trademark of Amersham Biosciences Corp. Cy™ dyes are subject to proprietary rights of Amersham Biosciences Corp and Carnegie Mellon University and are made and sold under license from Amersham Biosciences Corp only for research and in vitro diagnostic use. Any other use requires a commercial sublicense from Amersham Biosciences Corp. 800 Centennial Avenue, Piscataway, NJ 08855-1327, USA.

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

BD, BD Logo and all other trademarks are property of Becton, Dickinson and Company. © 2010 BD  
23-11911-00

## BD Biosciences

2350 Qume Drive  
San Jose, CA 95131  
US Orders: 877.232.8995  
answers@bd.com  
bdbiosciences.com