

Mouse IL-21 ELISPOT

Catalog Number: 88-7210

Also known as: Interleukin-21

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

Product Information

Contents: Mouse IL-21 ELISPOT
Catalog Number: 88-7210



Temperature Limitation: Store at 2-8°C.

Batch Code: Refer to vial

Use By: Refer to vial

Description

IL-21 is a 17 kDa immunomodulatory cytokine produced mainly by NKT, T helper (Th) 17 and T follicular helper (TFH) cells. In TFH cells, IL-21 expression leads to autocrine signaling through the IL-21 receptor (IL-21R) and STAT3, which leads to additional transcriptional activation by Bcl-6. As with IFN gamma for Th1, IL-4 for Th2 cells, and IL-17A for Th17 cells, IL-21 is critical for TFH effector function. This cytokine plays a role in T cell-dependent B cell differentiation into plasma cells and memory cells, stimulation of IgG production and induction of apoptotic signaling in naïve B cells.

In Th17 cells, IL-21 expression and autocrine feedback through STAT3, IRF4 and ROR gamma t lead to upregulation of the IL-23R, thereby preparing Th17 cells for maturation and maintenance by the inflammatory cytokine IL-23. While upregulating IRF4 and ROR gamma t, IL-21 also mediates the downregulation of Foxp3. High levels of IL-21 are present in chemically-induced colitis models. IL-21-deficient mice are protected from developing colitis upon chemical treatment by their inability to upregulate Th17-associated molecules.

Components

Capture Antibody. Pre-titrated, Functional Grade (low endotoxin) purified antibody

Detection Antibody. Pre-titrated, biotin-conjugated antibody

ELISA/ELISPOT Coating Buffer. This Ready-Set-Go! ELISPOT Set may contain ELISA/ELISPOT Coating Buffer Powder (Reconstitute to 1L with dH2O and filter (0.22 µM)) or 10X PBS ELISPOT Coating Buffer (Dilute 1 part 10X Buffer into 9 parts dH2O and filter with 0.22 µM).

Assay Diluent. 5X Concentrated

Detection Enzyme. Pre-titrated Avidin-HRP

Certificate of Analysis. Lot-specific instructions for the dilution of antibodies and enzyme

Applications Reported

This ELISPOT set is for the high resolution frequency analysis of IL-21 secreting cells.

Applications Tested

This mouse IL-21 ELISPOT Ready-SET-Go! set contains all of the necessary reagents for performing enzyme-linked immunosorbent spot (ELISPOT) assays for high resolution frequency analysis of IL-21 secreting cells. The reagents in this set have been pre-titrated for optimal spot development. Millipore Multiscreen HTS 96-well filtration plates are recommended, but not included, for use in this assay.

This assay has been validated for the detection of endogenous IL-21 using Th17-polarized Balb/c splenocytes. Splenocytes were cultured in the presence of Anti-Mouse CD3e Functional Grade Purified (cat. 16-0031), Anti-Mouse CD28 Functional Grade Purified (cat. 16-0281), Anti-Mouse IL-2 Functional Grade Purified (cat. 16-7022), Anti-Mouse IL-4 Functional Grade Purified (cat. 16-7041), Anti-Mouse IFN gamma Functional Grade Purified (cat. 16-7311), Human TGF beta 1 Recombinant Protein (cat. 14-8348), and Mouse IL-6 Recombinant Protein (cat. 14-8061) for three days to induce Th17 polarization. The cells were then stimulated with PMA and Ionomycin and transferred to coated plates for the analysis of IL-21 secretion.

References

Konforte D, Paige CJ. Interleukin-21 regulates expression of the immediate-early lytic cycle genes and proteins in Epstein-Barr Virus infected B cells. *Virus Res.* 2009 Sep;144(1-2):339-43.

Not for further distribution without written consent.

Copyright © 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

Tel: 888.999.1371 or 858.642.2058 • Fax: 858.642.2046 • thermofisher.com/ebioscience • info@ebioscience.com

Mouse IL-21 ELISPOT

Catalog Number: 88-7210

Also known as: Interleukin-21

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

Pot C, Jin H, Awasthi A, Liu SM, Lai CY, Madan R, Sharpe AH, Karp CL, Miaw SC, Ho IC, Kuchroo VK. Cutting edge: IL-27 induces the transcription factor c-Maf, cytokine IL-21, and the costimulatory receptor ICOS that coordinately act together to promote differentiation of IL-10-producing Tr1 cells. *J Immunol.* 2009 Jul 15;183(2):797-801.

Elsaesser H, Sauer K, Brooks DG. IL-21 is required to control chronic viral infection. *Science.* 2009 Jun 19;324(5934):1569-72.

Bauquet AT, Jin H, Paterson AM, Mitsdoerffer M, Ho IC, Sharpe AH, Kuchroo VK. The costimulatory molecule ICOS regulates the expression of c-Maf and IL-21 in the development of follicular T helper cells and TH-17 cells. *Nat Immunol.* 2009 Feb;10(2):167-75.

Nurieva R, Yang XO, Martinez G, Zhang Y, Panopoulos AD, Ma L, Schluns K, Tian Q, Watowich SS, Jetten AM, Dong C. Essential autocrine regulation by IL-21 in the generation of inflammatory T cells. *Nature.* 2007 Jul 26;448(7152):480-3.

Related Products

12-7211 eBioscience™ Anti-Mouse IL-21 PE (FFA21)

16-7211 eBioscience™ Anti-Mouse IL-21 Functional Grade Purified (FFA21)

88-8210 Mouse IL-21 Uncoated ELISA Kit

Not for further distribution without written consent.

Copyright © 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

Tel: 888.999.1371 or 858.642.2058 • Fax: 858.642.2046 • thermofisher.com/ebioscience • info@ebioscience.com

ELISPOT Protocol

Protocol: ELISPOT

Materials Provided

- Please refer to the Certificate of Analysis (C of A) for components

Other Materials Needed (see note at the end of the protocol)

- 96-Well PVDF Membrane ELISPOT Plates (Millipore, Cat. No. MAIPS4510)
- AEC (3-amino-9-ethyl carbazole) Substrate (Sigma, Cat. No. A-5754)
- Distilled water (dH₂O)
- ELISPOT Wash Buffer: 1X PBS, with 0.05% Tween-20 (or Thermo Fisher ELISA/ELISPOT Wash Buffer Powder, Cat. No. 00-0400)
- Complete RPMI-1640
- 1X PBS

Instruments

- Pipettes and pipettors
- Refrigerator
- Incubator
- Laminar Flow Hood
- Plate Washer: Wash bottle or automated wash machine
- ELISPOT plate reader or dissecting microscope for visual inspection

Time Requirements

- 1 overnight incubation
- 1-2 day cell activation
- 3-5 hour washing, antibody incubations and color development

Experimental Procedure

Aseptic Steps:

Note: Use sterile buffers and aseptic technique; perform all steps in a laminar flow hood.

1. Dilute Functional Grade purified capture antibody in sterile ELISA/ELISPOT Coating Buffer, as noted on Certificate of Analysis which is included with the reagent set. Coat ELISPOT plate with 100 μ L/well of capture antibody solution. Incubate at 2-8°C overnight.
2. Decant or aspirate coating antibody from plate.
3. Wash plates 2 times with 200 μ L/well of sterile ELISA/ELISPOT Coating Buffer. Decant.
4. Block plate with 200 μ L/well of complete RPMI-1640 at room temperature for 1 hour. Decant or aspirate plate.
5. Aliquot mitogen, antigen, or controls diluted in complete RPMI-1640 to appropriate wells at 100 μ L/well. Aliquot cells at desired densities (e.g., 1×10^5 /mL - 2×10^6 /mL) at 100 μ L/well and incubate at 37°C, in a 5% CO₂ humidified incubator for 24-48 hours.

Note: Optimal kinetics and cell densities vary with target cytokine, treatment, and cell type and must be empirically determined. See references. Cells can be diluted in a sterile tissue culture plate starting at 2×10^6 /well in triplicate wells with a series of 1:3 or 1:4 serial dilutions down the plate, and then transferred to the ELISPOT plate.

Non-Aseptic Steps:

6. Decant cells and medium from plates. Wash plate 3 times with ELISA/ELISPOT Wash Buffer.
7. Dilute biotinylated detection antibody in Assay Diluent according to instructions on the Certificate of Analysis provided with the reagent set. Add 100 μ L/well of detection antibody solution. Incubate at room temperature for 2 hr (or at 2-8°C overnight).

8. Decant antibody solution. Wash 4 times with ELISA/ELISPOT Wash Buffer. Allow wells to soak for 1 minute for each wash.
9. Dilute Avidin-HRP reagent in Assay Diluent according to instructions on the Certificate of Analysis provided with the reagent set. Add 100 μ L/well of Avidin-HRP solution and incubate at room temperature for 45 minutes.
10. Decant Avidin-HRP solution. Wash plate 3 times with ELISA/ELISPOT Wash Buffer, and then 2 times with 1X PBS (no Tween-20).
11. Add 100 μ L/well of freshly-prepared AEC Substrate Solution and develop at room temperature for 10-60 minutes; monitor development of spots.
12. Stop the substrate reaction by washing wells 3 times with 200 μ L/well of distilled water.
13. Air-dry the plate. Count spots using a dissecting microscope or automated ELISPOT plate reader. Store plates in the dark prior to reading.

Solutions

ELISA/ELISPOT Coating Buffer Powder:

- Reconstitute powder to 1 L in dH₂O; filter sterilize using a 0.22 μ M filter

Complete RPMI-1640:

- RPMI-1640 with 10% Fetal Bovine Serum and 1% Penicillin/Streptomycin/L-Glutamine

Assay Diluent (supplied as 5X)

- Dilute 5X solution to 1X in dH₂O

ELISA/ELISPOT Wash Buffer:

- 1X PBS with 0.05% Tween-20 (0.5 mL Tween-20 in 1 L PBS) or Thermo Fisher ELISA Wash Buffer Powder (Cat. No. 00-0400)

AEC (3-amino-9-ethyl carbazole) Substrate Solution:

- AEC Stock Solution: Dissolve 100 mg of AEC in 10 mL of N,N Dimethylformamide (DMF; Pierce, Cat. No. 20672)
- Add 333 μ L of AEC Stock Solution to 10 mL of 0.1 M Acetate Solution (pH 5.0) (see below for recipe). Filter through a 0.45 μ m filter.
- Just before use, add 5 μ L of 30% H₂O₂. Mix and use immediately.

0.1 M Acetate Solution (pH 5.0):

- Combine 148 mL of 0.2 M acetic acid (11.55 mL glacial acetic acid per 1 L of dH₂O) with 352 mL of 0.2 M sodium acetate (27.2 g sodium acetate per 1 L of dH₂O).
- QS to 1 L with dH₂O. Adjust pH to 5.0.

Documentation and support

Customer and technical support

Visit thermofisher.com/support for the latest in services and support, including:

- Worldwide contact telephone numbers
- Product support, including:
 - Product FAQs
 - Software, patches, and updates
- Order and web support
- Product documentation, including:
 - User guides, manuals, and protocols
 - Certificates of Analysis
 - Safety Data Sheets (SDSs; also known as MSDSs)

Note: For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

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.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at thermofisher.com/support.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

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.

Corporate entity: Life Technologies | Carlsbad, CA 92008 USA | Toll Free in USA 1.800.955.6288

©2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. All other trademarks are properties of their respective owners.

For support visit thermofisher.com/support or email techsupport@lifetech.com

thermofisher.com

23 January 2017

ThermoFisher
SCIENTIFIC