


## Mouse IL-4 ELISPOT

**Catalog Number:** 88-7844

**Also known as:** Interleukin-4

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

### Product Information

 **Contents:** Mouse IL-4 ELISPOT  
**Catalog Number:** 88-7844



**Temperature Limitation:** Store at 2-8°C.



**Batch Code:** Refer to vial



**Use By:** Refer to vial



**Contains sodium azide**

### Description

This Mouse IL-4 ELISPOT Ready-SET-Go! reagent set contains the necessary reagents for performing enzyme linked immunosorbent spot (ELISPOT) assays for high resolution frequency analysis of IL-4-secreting cells. This ELISPOT reagent set is pre-titrated for optimal spot development.

### Components

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

**Detection Antibody.** Pre-titrated, biotin-conjugated antibody

**ELISA/ELISPOT Coating Buffer Powder.** 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 dilution of antibodies and enzyme

### References

Anis MM, Fulton SA, et al. 2007. Modulation of naive CD4+ T-cell responses to an airway antigen during pulmonary mycobacterial infection. Infect Immun. 75(5):2260-8. (ELISPOT kit, PubMed)

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](http://thermofisher.com/ebioscience) • [info@ebioscience.com](mailto: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 (dH<sub>2</sub>O)
- 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  $\mu$ 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  $\mu$ L/well of sterile ELISA/ELISPOT Coating Buffer. Decant.
4. Block plate with 200  $\mu$ 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  $\mu$ L/well. Aliquot cells at desired densities (e.g.,  $1 \times 10^5$ /mL -  $2 \times 10^6$ /mL) at 100  $\mu$ L/well and incubate at 37°C, in a 5% CO<sub>2</sub> 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  $2 \times 10^6$ /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  $\mu$ 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  $\mu$ 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  $\mu$ 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  $\mu$ 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 dH<sub>2</sub>O; filter sterilize using a 0.22  $\mu$ 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 dH<sub>2</sub>O

### 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  $\mu$ 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  $\mu$ m filter.
- Just before use, add 5  $\mu$ L of 30% H<sub>2</sub>O<sub>2</sub>. 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 dH<sub>2</sub>O) with 352 mL of 0.2 M sodium acetate (27.2 g sodium acetate per 1 L of dH<sub>2</sub>O).
- QS to 1 L with dH<sub>2</sub>O. Adjust pH to 5.0.

## Documentation and support

### Customer and technical support

Visit **[thermofisher.com/support](http://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](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html)**. If you have any questions, please contact Life Technologies at **[thermofisher.com/support](http://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](http://thermofisher.com/support)** or email **[techsupport@lifetech.com](mailto:techsupport@lifetech.com)**

**[thermofisher.com](http://thermofisher.com)**

23 January 2017

**ThermoFisher**  
SCIENTIFIC