

---

## Batf3 Polyclonal Antibody, FITC Conjugated

(Catalog # A50087)

---

### Background

AP-1 family transcription factor that controls the differentiation of CD8+ thymic conventional dendritic cells in the immune system. Acts via the formation of a heterodimer with JUN family proteins that recognizes and binds DNA sequence 5'-TGA[CG]TCA-3' and regulates expression of target genes. Required for development of CD8-alpha+ classical dendritic cells (cDCs) and related CD103+ dendritic cells that cross-present antigens to CD8 T-cells and produce interleukin-12 (IL12) in response to pathogens.

### Description

Batf3 Polyclonal Antibody, FITC Conjugated. FITC. Raised in: Rabbit.

### Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

### Specificity

Rat

### Isotype

IgG

### Uniprot ID

P97876

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Rat Basic leucine zipper transcriptional factor ATF-like 3 protein (1-133AA)

### Storage

Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

B-ATF-3, Jun dimerization protein 1, JDP-1, Jdp1, Batf3

### Application

ELISA; Recommended dilution: ELISA 1:100-1:500

---