



Anti-Mouse CD3 ϵ (Biotinylated)

CATALOG NO.: 54726

CLONE: 145-2C11

BACKGROUND:

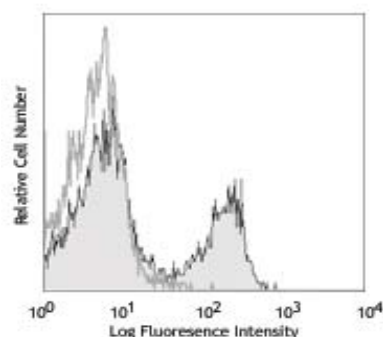
CD3 ϵ is a 20 kD transmembrane protein, also known as CD3 or T3. It is a member of the Ig superfamily and primarily expressed on T cells, NK-T cells, and at different levels on thymocytes during T cell differentiation. The 145-2C11 antibody is useful for *in vitro* blocking and activation assays, as well as apoptosis induction and *in vivo* T cell depletions.

APPLICATION:

Applications include flow cytometry, immunohistochemistry¹⁴, immunofluorescence microscopy⁵, western blotting⁴, immunoprecipitation¹, immunohistochemical staining¹⁴ of acetone-fixed frozen sections and zincfixed paraffin-embedded sections, Western blotting⁴, complement-mediated cytotoxicity⁶, *in vitro* and *in vivo* stimulation of T cells^{1,2,7,12}, immunofluorescent staining⁵, and *in vivo* T cell depletions⁸⁻¹⁰. Recommended usage: Each lot of this antibody is quality control tested by immunofluorescent staining with flow cytometric analysis. For immunofluorescent staining, the suggested use of this reagent is $\leq 0.25 \mu\text{g}$ per 10^6 cells in 100 μl . It is recommended that the reagent be titrated for optimal performance for each application.

REFERENCES:

1. Leo, O., *et al.*, 1987. *PNAS* 84:1374.
2. Kruisbeek, A.M., *et al.*, 1991. *In Current Protocols in Immunology*. pp. 3.12.1 - 3.12.14.
3. Duke, R.C., *et al.*, 1995. *Current Protocols in Immunology*. pp. 3.17.1 - 3.17.33.
4. Salvadori, S., *et al.*, 1994. *J. Immunol.* 153:5176. 5. Payer, E., *et al.*, 1991. *J. Immunol.* 146:2536.
6. Jacobs, H., *et al.*, 1994. *Eur. J. Immunol.* 24:934.
7. Vossen, A.C.T.M., *et al.*, 1995. *Eur. J. Immunol.* 25:1492.
8. Henrickson, M., *et al.*, 1995. *Transplantation* 60:828.
9. Kinnaert, P., *et al.*, 1996. *Transpl. Int.* 9:386.
10. Han, W.R., *et al.*, 1999. *Transpl. Immunol.* 7:207.
11. Miescher, G.C., *et al.*, 1989. *Immunol. Lett.* 23:113.
12. Terrazas, L. I., *et al.*, 2005. *Intl. J. Parasitology*. 35:1349.
13. Ko, S-Y., *et al.*, 2005. *J. Immunol.* 175:3309.
14. Podd, B.S., *et al.*, 2006. *J. Immunol.* 176: 6532.



C57BL/6 mouse splenocytes stained with purified 145-2C11, followed by anti-Armenian hamster IgG FITC

This product is for in vitro research purposes only.

STORAGE:

This monoclonal antibody was purified by affinity chromatography. The vial contains 0.25 ml antibody (at 200 $\mu\text{g/ml}$) in 1X PBS (pH 7.2) with 0.09% sodium azide. Stored at 4°C. for up to 1 year. Avoid repeated freezing and thawing.

RELATED PRODUCTS:

- Streptavidin; HiLyte Fluor 680 labeled, Cat # **60659-H680**
- Streptavidin; HiLyte Fluor 750 labeled, Cat # **60659-H750**
- Streptavidin; HiLyte Fluor 488 labeled, Cat # **60665**
- Streptavidin; HiLyte Fluor 647 labeled, Cat # **60667**
- Streptavidin; HiLyte Fluor 555 labeled, Cat # **60666**
- Streptavidin; HiLyte Fluor TR labeled, Cat # **60671**
- Streptavidin; HiLyte Plus 555 labeled, Cat # **60672-Plus555**
- Streptavidin; HiLyte Plus 647 labeled, Cat # **60672-Plus647**
- Streptavidin; HiLyte Plus 750 labeled, Cat # **60672-Plus750**