

## PRODUCT INSERT

# MOUSE anti-MOUSE CD72.2

Product Code	Form	Volume	Antibody*	Excitation (nm)	Peak Emission (nm)	Matching Isotype Controls	
MM5800	Purified	1.0 ml	200 µg	N/A	N/A	Mouse IgM Purified	Code MGM00
MM5815	Biotin	1.0 ml	100 µg	N/A	N/A	Mouse IgM Biotin	Code MGM15
MM5815-3	Biotin	3.0 ml	300 µg				

## PRODUCT DESCRIPTION

Mouse monoclonal antibody to mouse CD72.2

**Clone:** CT-72.2

**Isotype:** Mouse IgM

**Lot No.:** See label      **Expiration:** See label

**Buffer:** Phosphate buffered saline (PBS)

**Preservatives:** 0.1% *sodium azide*. Sodium azide is an extremely toxic and dangerous compound particularly when combined with acids or metals. Solutions containing sodium azide should be disposed of properly.

**Stabilizer:** For conjugated products only, a highly purified grade of BSA has been added as a stabilizing protein..

## STORAGE AND HANDLING

Store reagents at 2-8°C. Light exposure should be avoided with fluorochrome conjugated reagents. Use dim light during handling, incubation with cells and prior to analysis. If the reagent is being diluted, it is recommended that only the quantity to be used within one week be diluted.

Using formalin to fix cells following immunofluorescent staining may cause the degradation of tandem fluorochromes. Cells stained with TRI-COLOR®, PE-Cy7, PE-TR or APC-Cy7 should be analyzed by flow cytometry within 18 hours following fixation.

## PRODUCT CHARACTERIZATION

**Antigen Specificity:** The CT-72.2 monoclonal antibody (mAb) reacts with the CD72 alloantigen CD72.2, a B-cell surface protein that is encoded by the *Cd72<sup>b</sup>* allele. CD72.2 is expressed on B lymphocytes and subsets of T cells. Mouse strains expressing CD72.2 include BALB/c, CBA/ca, CBA/N, C3H/He, C57BL/-, 129/- NZB/- and PL/J. The CT-72.2 mAb can be used in immunostaining for flow cytometry.

## PRODUCT QUALITY CONTROL

Every lot is tested by flow cytometry using freshly harvested mouse splenocytes. From this testing it is recommended that between 0.1 and 0.25µg of antibody be used per 1 x 10<sup>6</sup> cells in a 100µl staining volume. Because conditions may vary, it is recommended that each investigator determine the optimal amount of antibody to be used for their application.

## REFERENCES

1. Ying, H., J. I. Healy, C. C. Goodnow, and J. R. Parnes. 1998. Regulation of mouse CD72 gene expression during B lymphocyte development. *J. Immunol.* 161: 4760-4767.

\* The amount of antibody is determined by measuring the optical density using a spectrophotometer. The antibody titer is verified by immunofluorescent staining and flow cytometric analysis.

**FOR RESEARCH USE ONLY. . . NOT FOR THERAPEUTIC OR IN VITRO DIAGNOSTIC USE**