

ANTI-HUMAN CD89 PE

**PE-Conjugated Mouse Monoclonal Antibody Against Human CD89
Clone A59**

Catalog # 10530

100 tests

SPECIFICITY:

CD89 is the Fc receptor for immunoglobulin A (IgA, Fc α R) and is expressed on promyelocytes, promonocytes, mature neutrophils, monocytes/ macrophages and on activated eosinophils. Anti-CD89 clone A59 does not block the binding of IgA to the Fc α R.

CLONE: A59

ISOTYPE: IgG₁ κ (mouse)

FORMAT:

PE-conjugated antibody in 2 mL aqueous buffered solution containing bovine serum albumin and 0.09% (w/v) sodium azide.

STABILITY AND STORAGE:

Store at 4°C. Do not freeze. Product is stable for at least 6 months.

APPLICATIONS AND DIRECTIONS FOR USE:

Flow cytometry:

Recommended amount per 1x10⁶ cells in a volume of 100 μ L: 20 μ L
Appropriate conditions should be established for each application.

Cell separation:

Positive selection of CD89⁺ cells with StemCell's reagents for immunomagnetic cell separation. Please contact us for more information.

**NOT FOR CLINICAL USE; INCLUDING IN VITRO DIAGNOSTIC USE, AND EX VIVO OR IN VIVO THERAPEUTIC USE IN CLINICAL TRIALS OR IN CLINICAL PRACTISE.
THIS REAGENT IS FOR RESEARCH ONLY.**

Hazardous Ingredient: Sodium Azide. Avoid exposure to skin and eyes, ingestion and contact with heat, acids and metals. Wash exposed skin with soap and water. Flush eyes with water. Dilute with running water before discharging into plumbing.

REFERENCES:

1. Barclay AN, Brown MH, Law SKA, McKnight AJ, Tomlinson MG, van der Merwe PA, eds. 1997. The Leukocyte Antigens Facts Book, 2nd Edition, CD89. Academic Press, NY, p 351.
2. Kishimoto T et al. Eds.1998. Leukocyte Typing VI. White Cell Differentiation Antigens.. Garland Publishing Inc, New York NY, p1178.