

ANTI-HUMAN CD36

**Monoclonal Antibody Against Human CD36
Clone FA6-152**

Catalog # 01436

100 mg

PRODUCT DESCRIPTION:

CD36 is a 88 kDa glycoprotein receptor for extracellular matrix proteins such as collagen and thrombospondin. CD36 is expressed on monocyte, macrophages, platelets and erythrocyte precursors. CD36 is also expressed on endothelial and epithelial cells, adipocytes and activated keratinocytes.

CLONE: FA6-152

ISOTYPE: IgG₁, κ (mouse)

FORMULATION:

100 µg in 0.1 mL (1 mg/mL) of Phosphate-buffered Saline (PBS). Does not contain sodium azide or albumin. The antibody was purified from tissue hybridoma culture supernatant by Protein G affinity chromatography.

It should be kept in mind that this product is a biological reagent, and as such cannot be completely characterized or quantified. Some variability is unavoidable.

STABILITY AND STORAGE:

Store at 4°C. Do not freeze. Product is stable for at least 2 years.

Contents guaranteed sterile if seal is not tampered with.

APPLICATIONS AND DIRECTIONS FOR USE:

Flow cytometry:

Recommended amount per 1×10^6 cells in a volume of 100 µL: ≤ 1 µg (1 µL)

Cell separation:

Positive selection or depletion of human CD36⁺ cells with StemCell's reagents for immunomagnetic cell separation. Please contact us for more information.

**THIS REAGENT IS FOR RESEARCH USE ONLY.
IT IS NOT TO BE ADMINISTERED TO HUMANS.**

REFERENCES:

1. CD36. In: The Leukocyte Antigen FactsBook. Barclay AN, Brown MH, Law SKA et al Eds. Academic Press Inc., San Diego, USA, 1997, pp221-223.
2. CD36. In: Leukocyte Typing VII. Mason D et al. Eds., Oxford University Press, Oxford UK, 2002, pp781-783