

**APC/iFluor™ 750 Anti-human CD53
Antibody *HI36***Catalog number: 105311G0, 105311G1, 105311G2
Unit size: 25 tests, 100 tests, 500 tests**Product Details**

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	0.1 mg/mL
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG3
Immunogen	CD53 (Tetraspanin-25, MOX44)
Clone	HI36
Conjugate	APC/iFluor™ 750

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with APC/iFluor™ 750 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	APC/iFluor™ 750
Excitation Wavelength	754 nm
Emission Wavelength	793 nm

Applications

The HI36 monoclonal antibody binds to human CD53, a 35 - 42 kD member of the tetraspanin/transmembrane-4 superfamily typically expressed on the surface of leukocytes, B cells, dendritic cells, osteoblasts and thymocytes. In many organisms, CD53 plays a role in the upregulation of myoblast fusion, and is associated with a variety of biologically interesting macromolecules/ligands, for instance, VLA-4 and HLA-DR. CD53 is a relatively rare antibody target, with fewer than 600 publications in the last decade. Even still, CD53 is vital to costimulatory molecules research, commonly serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through

affinity chromatography and conjugated to APC/iFluor™ 750 (ex/em = 754/793 nm).