

iFluor™ 488 Anti-human CD243 Antibody
UIC2Catalog number: 12430050, 12430051
Unit size: 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 IgG2a kappa
Immunogen	CD243 (MDR-1)
Clone	UIC2
Conjugate	iFluor™ 488

Biological Properties

Appearance	Orange-red liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 488 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 488
Excitation Wavelength	491 nm
Emission Wavelength	516 nm

Applications

The UIC2 monoclonal antibody recognizes human CD243, a 170 kD member of the Member of ABC transporter family typically located on the surface of hematopoietic stem cells, T cells, natural killer cells and B cells. In some organisms, CD243 is involved in the positive regulation of anion channel activity, and is associated with a variety of biologically interesting macromolecules/ligands. CD243 is a relatively rare antibody

target, with fewer than 100 publications in the last decade. Even still, CD243 is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, specifically in the study of immunology, cancer biomarkers and cell biology. This antibody was purified through affinity chromatography and conjugated to iFluor™ 488 (ex/em = 491/516 nm). It is compatible with the 488 nm laser and 530/30 nm bandpass filter (for example, as in the BD FACSCelesta™).