

**mFluor™ Blue 570 Anti-human CD243
Antibody *UIC2***Catalog number: 124300T0, 124300T1
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	mFluor™ Blue 570

Biological Properties

Appearance	Red liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ Blue 570 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	mFluor™ Blue 570
Excitation Wavelength	553 nm
Emission Wavelength	565 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 mFluor™ Blue 570 (ex/em = 553/565 nm). It is compatible with the 561 nm laser and 583/24 nm bandpass filter (for example, as in the Luminex Amnis CellStream).