

**mFluor™ Red 700 Anti-human CD158d
Antibody *mAb#33***Catalog number: 115810V0, 115810V1
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 IgG1 kappa
Immunogen	CD158d (KIR2DL4)
Clone	mAb#33
Conjugate	mFluor™ Red 700

Biological Properties

Appearance	Dark blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ Red 700 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	mFluor™ Red 700
Excitation Wavelength	680 nm
Emission Wavelength	695 nm

Applications

The mAb#33 monoclonal antibody binds to human CD158d, a 45 - 50 kD transmembrane glycoprotein frequently expressed on the surface of natural killer cells. In certain organisms, CD158d negatively regulates natural killer cell mediated cytotoxicity, is an enhancer of cellular senescence and positively regulates natural killer cell cytokine production. From a research standpoint, it is of biological interest due to its

association with essential macromolecules/ligands like HLA-Bw4. CD158d is a relatively rare antibody target, with fewer than 200 publications in the last decade. Even still, CD158d has been widely used in immunology and innate immunity research, often serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to mFluor™ Red 700 (ex/em = 680/695 nm). It is compatible with the 642 nm laser and 664/20 nm bandpass filter (for example, as in the Luminex Guava easyCyte).