

iFluor™ 555 Anti-human CD178 Antibody
NOK-1Catalog number: 11780090, 11780091
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
Immunogen	CD178 (Fas ligand, CD95L, FASL, TNFSF6)
Clone	NOK-1
Conjugate	iFluor™ 555

Biological Properties

Appearance	Red liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 555 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 555
Excitation Wavelength	557 nm
Emission Wavelength	570 nm

Applications

The NOK-1 monoclonal antibody binds to human CD178, a 38 - 42 kD member of the TNF superfamily commonly found on the surface of dendritic cells, natural killer cells, testis, neutrophils and T cells. CD178 is associated with a variety of biologically interesting macromolecules/ligands, for example, DcR3. CD178 is a relatively rare antibody target, with fewer than 500 publications in the last decade. Even

still, CD178 is often used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of cell biology, apoptosis/tumor suppressors/cell death and neuroscience. This antibody was purified through affinity chromatography and conjugated to iFluor™ 555 (ex/em = 557/570 nm). It is compatible with the 561 nm laser and 585/20 nm bandpass filter (for example, as in the BD Special Order LSRFortessa™ Cell Analyzer).