

XFD680 Anti-human CD178 Antibody *NOK-1*

Catalog Number: 11780180, 11780181

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	Lot specific (please consult certificate of analysis for given lot)
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	AF680

Biological Properties

Appearance	Blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF680 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging
Recommended Dilutions	For flow cytometry applications, the suggested concentration is at 5 uL/million cells in 100 uL staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined. <i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i>

Spectral Properties

Conjugate AF680

Excitation Wavelength 681 nm

Emission Wavelength 704 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 XFD680 (ex/em = 681/704 nm). XFD680 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 680 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 642 nm laser and 702/85 nm bandpass filter (for example, as in the Luminex Amnis FlowSight).