

**PE/XFD700 Anti-human CD95 Antibody \*EOS9.1\***

Catalog Number: 109501Q0,  
109501Q1, 109501Q2  
Unit Size: 25 tests, 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 IgM kappa
Immunogen	CD95 (Fas, TNFRSF6, APO-1, FASLG)
Clone	EOS9.1
Conjugate	PE/AF700

**Biological Properties**

---

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/AF700 under optimal conditions
Application	Flow Cytometry (FACS)
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	PE/AF700
-----------	----------

Excitation Wavelength 565 nm

Emission Wavelength 721 nm

## Applications

---

EOS9.1 is an anti-human monoclonal antibody that forms an immune complex with the CD95 antigen. CD95 (alternatively called APO-1 or TNFRSF6) is a 45 kD member of the TNFR superfamily that is found on the surface of cells like NK cells, T cells and granulocytes. CD95 is associated with a variety of biologically interesting macromolecules/ligands, for instance, CD178 (fas ligand). CD95 is a moderately popular antibody target, with over 12000 publications in the last decade. CD95 is commonly used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of cell biology. This antibody was purified through affinity chromatography and conjugated to PE/XFD700 (ex/em = 566/721 nm). XFD700 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 700 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 561 nm laser and 695/40 nm bandpass filter (for example, as in the Agilent Technologies NovoCytte Quanteon).