

iFluor™ 810 Anti-human CD95 Antibody
EOS9.1Catalog number: 10950000, 10950001
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 IgM kappa
Immunogen	CD95 (Fas, TNFRSF6, APO-1, FASLG)
Clone	EOS9.1
Conjugate	iFluor™ 810

Biological Properties

Appearance	Green liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 810 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 810
Excitation Wavelength	811 nm
Emission Wavelength	822 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 iFluor™ 810 (ex/em = 811/822 nm).