

**mFluor™ Blue 570 Anti-human CD95  
Antibody \*EOS9.1\***Catalog number: 109500T0, 109500T1  
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          | mFluor™ Blue 570                  |

**Biological Properties**

|             |                                                                                                                 |
|-------------|-----------------------------------------------------------------------------------------------------------------|
| Appearance  | Red liquid                                                                                                      |
| Preparation | Antibody purified by affinity chromatography and then conjugated with mFluor™ Blue 570 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                                     |

**Spectral Properties**

|                       |                  |
|-----------------------|------------------|
| Conjugate             | mFluor™ Blue 570 |
| Excitation Wavelength | 553 nm           |
| Emission Wavelength   | 565 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 mFluor™ Blue 570 (ex/em = 553/565 nm). It is compatible with the 561 nm laser and 577/35 nm bandpass filter (for example, as in the Luminex Amnis ImageStream).