

**TRITC Anti-human CD178 Antibody \*NOK-1\***Catalog number: 117801H0, 117801H1  
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          | TRITC                                   |

**Biological Properties**

|             |  |
|-------------|--|
| Preparation | Antibody purified by affinity chromatography and then conjugated with TRITC under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging  |

**Spectral Properties**

|                       |        |
|-----------------------|--------|
| Conjugate             | TRITC  |
| Excitation Wavelength | 544 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 TRITC (ex/em = 544/570 nm). It is compatible with the 532 nm laser and 583/24 nm bandpass filter (for example, as in the Luminex Amnis

CellStream).