

**PE/Texas Red® Anti-human/ dog CD132  
Antibody \*TUGh4\***Catalog number: 113201R0, 113201R1, 113201R2  
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      | 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, dog                    |
| Class              | Primary                       |
| Clonality          | Monoclonal                    |
| Host               | Rat                           |
| Isotype            | Rat IgG2b kappa               |
| Immunogen          | CD132 (Common $\gamma$ chain) |
| Clone              | TUGh4                         |
| Conjugate          | PE/Texas Red®                 |

**Biological Properties**

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

**Spectral Properties**

|                       |               |
|-----------------------|---------------|
| Conjugate             | PE/Texas Red® |
| Excitation Wavelength | 567 nm        |
| Emission Wavelength   | 615 nm        |

**Applications**

The TUGh4 monoclonal antibody binds to human/ dog CD132, a 64 - 70 kD member of the Ig superfamily frequently located on the surface of macrophages and B cells. CD132 is a component of important cellular pathways, namely, the interleukin-21-mediated signaling pathway, interleukin-9-mediated signaling pathway and interleukin-2-mediated signaling pathway. In addition, in certain organisms, it is an enhancer of phagocytosis. From a research standpoint, it is of biological interest due to its association with important macromolecules/ligands such as CD124, JAK1, CD25 and lck. CD132 is a relatively rare antibody target, with fewer than 800 publications in the last decade. Even still, CD132 has

a variety of applications in immunology research, frequently serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to PE/Texas Red® (ex/em = 567/615 nm). It is compatible with the 561 nm laser and 615/20 nm bandpass filter (for example, as in the Agilent Technologies NovoCyt<sup>e</sup> Quanteon).