

**PerCP/Cy5.5 Anti-human/ dog CD132
Antibody *TUGh4***Catalog number: 113201T0, 113201T1, 113201T2
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 γ chain)
Clone	TUGh4
Conjugate	PerCP/Cy5.5

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PerCP/Cy5.5 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PerCP/Cy5.5
Excitation Wavelength	489 nm
Emission Wavelength	679 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 PerCP/Cy5.5 (ex/em = 489/679 nm). It is compatible with the 488 nm laser and 670/30 nm bandpass filter (for example, as in the BD LSRFortessa™ Cell Analyzer).