

**FITC Anti-human/ dog CD132 Antibody
*TUGh4***Catalog number: 113201G0, 113201G1
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, dog
Class	Primary
Clonality	Monoclonal
Host	Rat
Isotype	Rat IgG2b kappa
Immunogen	CD132 (Common γ chain)
Clone	TUGh4
Conjugate	FITC

Biological Properties

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

Spectral Properties

Conjugate	FITC
Excitation Wavelength	491 nm
Emission Wavelength	516 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 Ick. 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 FITC (ex/em = 491/516 nm). It is compatible with the 488 nm laser and 525/40 nm bandpass filter (for example, as in the Beckman Coulter DxFLEx).