

XFD594 Anti-human CD150 Antibody *SLAM.4*

Catalog Number: 11500160, 11500161

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	Lot specific (please consult certificate of analysis for given lot)
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	CD150 (SLAM, IPO-3)
Clone	SLAM.4
Conjugate	AF594

Biological Properties

Appearance	Purple liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF594 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging
Recommended Dilutions	For flow cytometry applications, the suggested concentration is at 5 uL/million cells in 100 uL staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined. <i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i>

Spectral Properties

Conjugate AF594
Excitation Wavelength 590 nm
Emission Wavelength 618 nm

Applications

The SLAMF4 monoclonal antibody binds with human CD150, a 70 - 95 kD glycoprotein commonly located on the surface of B cells, dendritic cells, endothelial cells, T cells and Tregs. CD150 plays a role in important cellular pathways, in particular, the negative regulation of CD40 signaling pathway. In addition, in certain organisms, it is a negative regulator of tumor necrosis factor production, suppresses interleukin-6 production and plays a role in the upregulation of JNK cascade. From a research standpoint, it is of biological interest due to its association with critical macromolecules/ligands such as tyrosine phosphatase CD45. CD150 is a fairly uncommon antibody target, with a little more than 2700 publications in the last decade. Even still, CD150 is often used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of immunology. This antibody was purified through affinity chromatography and conjugated to XFD594 (ex/em = 590/618 nm). XFD594 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 594 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 592 nm laser and 610/30 nm bandpass filter (for example, as in the Luminex Amnis ImageStream).