

XFD350 Anti-human CD8 Antibody *HIT8a*

Catalog Number: 10080140, 10080141

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	CD8a (T8, Leu2)
Clone	HIT8a
Conjugate	AF350

Biological Properties

Appearance	Off-white liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF350 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 AF350
Excitation Wavelength 343 nm
Emission Wavelength 441 nm

Applications

The HIT8a monoclonal antibody binds to human CD8a, a 32 - 34 kD member of the Ig superfamily often located on the surface of dendritic cells, t cells and thymocytes. CD8 is a component of important cellular pathways, namely, the transmembrane receptor protein tyrosine kinase signaling pathway and cell surface receptor signaling pathway. From a research standpoint, it is of biological interest due to its association with key macromolecules/ligands like Lck and MHCI. CD8 is a very popular antibody target, with over 125000 publications in the last decade. CD8a is frequently 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 XFD350 (ex/em = 343/441 nm). XFD350 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 350 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 355 nm laser and 450/50 nm bandpass filter (for example, as in the BD LSRFortessa™ Cell Analyzer).