

XFD647 Anti-human CD102 Antibody *CBR-IC2/2*

Catalog Number: 11020170, 11020171

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 IgG2a
Immunogen	CD102 (ICAM2)
Clone	CBR-IC2/2
Conjugate	AF647

Biological Properties

Appearance	Dark blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF647 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 AF647

Excitation Wavelength 650 nm

Emission Wavelength 671 nm

Applications

CBR-IC2/2 is an anti-human monoclonal antibody that targets the CD102 antigen. CD102 (alternatively called ICAM2) is a 55 kD transmembrane glycoprotein that is expressed on the surface of cells like T cells and macrophages. CD102 is associated with a variety of biologically interesting macromolecules/ligands, for instance, LFA-1, integrin $\alpha\beta 2$ and CD11b/CD18. CD102 is a relatively rare antibody target, with fewer than 200 publications in the last decade. Even still, CD102 is often used in flow cytometry applications as a phenotypic marker for differentiation of cell types, specifically in the study of immunology. This antibody was purified through affinity chromatography and conjugated to XFD647 (ex/em = 650/671 nm). XFD647 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 647 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 642 nm laser and 664/20 nm bandpass filter (for example, as in the Luminex Guava easyCyte).