

iFluor™ 840 Anti-human CD102 Antibody
CBR-IC2/2Catalog number: 110200Q0, 110200Q1
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
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG2a
Immunogen	CD102 (ICAM2)
Clone	CBR-IC2/2
Conjugate	iFluor™ 840

Biological Properties

Appearance	Dark brown liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 840 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 840
Excitation Wavelength	836 nm
Emission Wavelength	879 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 iFluor™ 840 (ex/em = 836/879 nm).