

## APC/XFD750 Anti-human/ non-human primates CD103 Antibody \*Ber-ACT8\*

Catalog Number: 110301D0,  
110301D1, 110301D2  
Unit Size: 25 tests, 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, non-human primates
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1 kappa
Immunogen	CD103 (Integrin alpha E, ITGAE)
Clone	Ber-ACT8
Conjugate	APC/AF750

### Biological Properties

---

Preparation	Antibody purified by affinity chromatography and then conjugated with APC/AF750 under optimal conditions
Application	Flow Cytometry (FACS)
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	APC/AF750
-----------	-----------

Excitation Wavelength 651 nm

Emission Wavelength 785 nm

## Applications

---

Ber-ACT8 is an anti-human/ non-human primates monoclonal antibody that forms an immune complex with the CD103 antigen. CD103 (sometimes referred to as Integrin alpha E) is a 150 kD member of the integrin family that is found on the surface of cells such as T cells. CD103 is associated with a variety of biologically interesting macromolecules/ligands, namely, integrin  $\beta$ 7 and E-Cadherin. CD103 is a fairly uncommon antibody target, with a little more than 7500 publications in the last decade. Even still, CD103 is often used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of neuroscience, immunology and synaptic biology. This antibody was purified through affinity chromatography and conjugated to APC/XFD750 (ex/em = 756/785 nm). XFD750 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 750 (Alexa Fluor® is the trademark of Thermo Fisher).