

**mFluor™ Violet 450 Anti-human/ non-human primates CD103 Antibody \*Ber-ACT8\***Catalog number: 110300Z0, 110300Z1  
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, non-human primates
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1 kappa
Immunogen	CD103 (Integrin alpha E, ITGAE)
Clone	Ber-ACT8
Conjugate	mFluor™ Violet 450

**Biological Properties**

Appearance	Light yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ Violet 450 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

**Spectral Properties**

Conjugate	mFluor™ Violet 450
Excitation Wavelength	406 nm
Emission Wavelength	445 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 mFluor™ Violet 450 (ex/em = 406/445 nm). It is compatible with the 405 nm laser and 457/45 nm bandpass filter (for example, as in the Luminex Amnis FlowSight).