

PE/Texas Red® Anti-human/ non-human primates CD103 Antibody *Ber-ACT8*Catalog number: 110301R0, 110301R1, 110301R2
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	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	PE/Texas Red®

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/Texas Red® under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/Texas Red®
Excitation Wavelength	567 nm
Emission Wavelength	615 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 PE/Texas Red® (ex/em = 567/615 nm). It is compatible with the 561 nm laser and 610/20 nm bandpass filter (for example, as in the BD Special Order LSRFortessa™ Cell Analyzer).