

**APC/iFluor™ 700 Anti-human/ non-human primates CD103 Antibody \*Ber-ACT8\***Catalog number: 110301E0, 110301E1, 110301E2  
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          | APC/iFluor™ 700                 |

**Biological Properties**

|             |                                                                                                                |
|-------------|----------------------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography and then conjugated with APC/iFluor™ 700 under optimal conditions |
| Application | Flow Cytometry (FACS)                                                                                          |

**Spectral Properties**

|                       |                 |
|-----------------------|-----------------|
| Conjugate             | APC/iFluor™ 700 |
| Excitation Wavelength | 685 nm          |
| Emission Wavelength   | 710 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/iFluor™ 700 (ex/em = 685/710 nm).