

PE/Texas Red® Anti-human/ non-human primates CD83 Antibody *HB15e*Catalog number: 108301U0, 108301U1, 108301U2
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	CD83 (HB15)
Clone	HB15e
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

HB15e is an anti-human/ non-human primates monoclonal antibody that is specific for the CD83 antigen. CD83 (alternatively called HB15) is a 43 kD member of the Ig superfamily that is located on the surface of cells like B cells and dendritic cells. CD83 is associated with a variety of biologically interesting macromolecules/ligands, for instance, . CD83 is a fairly uncommon antibody target, with a little more than 5100 publications in the last decade. Even still, CD83 has a variety of applications in immunology research, frequently serving as a phenotypic marker for differentiating cell types in flow cytometric applications. 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 620/15 nm bandpass filter (for example, as in the Thermo Fisher Attune NxT).