

**iFluor™ 350 Anti-non-human primates/
human CD35 Antibody *E11***Catalog number: 10350010, 10350011
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	Non-human primates, human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1
Immunogen	CD35 (C3b/C4b receptor, Complement receptor type 1)
Clone	E11
Conjugate	iFluor™ 350

Biological Properties

Appearance	Off-white liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 350 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 350
Excitation Wavelength	345 nm
Emission Wavelength	450 nm

Applications

E11 is an anti-non-human primates/ human monoclonal antibody that recognizes the CD35 antigen. CD35 (alternatively called Complement receptor type 1) is a single-pass type I membrane protein that is found on the surface of cells such as erythrocytes. CD35 plays a role in key cellular pathways, for example, the negative regulation of complement activation, alternative pathway and complement activation, classical

pathway. Moreover, in certain organisms, it represses serine-type endopeptidase activity, plays a role in the downregulation of complement activation, alternative pathway and is involved in the positive regulation of serine-type endopeptidase activity. From a research standpoint, it is of biological interest due to its association with important macromolecules/ligands such as ic3, c4b, c3b and ic4. CD35 is a fairly uncommon antibody target, with a little more than 1800 publications in the last decade. Even still, CD35 has been widely used in immunology, cell biology and neuroinflammation research, commonly serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to iFluor™ 350 (ex/em = 345/450 nm). It is compatible with the 355 nm laser and 447/60 nm bandpass filter (for example, as in the Bio-Rad ZE5 Cell Analyzer).