

**iFluor™ 488 Anti-non-human primates/
human CD35 Antibody *E11***Catalog number: 10350050, 10350051
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™ 488

Biological Properties

Appearance	Orange-red liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 488 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 488
Excitation Wavelength	491 nm
Emission Wavelength	516 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™ 488 (ex/em = 491/516 nm). It is compatible with the 488 nm laser and 530/30 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte Quanteon).