

**PE/iFluor™ 750 Anti-human CD2 Antibody**  
**\*HIT11\***Catalog number: 100201R0, 100201R1, 100201R2  
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
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1
Immunogen	CD2 (LFA-2, Erythrocyte R, T11)
Clone	HIT11
Conjugate	PE/iFluor™ 750

**Biological Properties**

---

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/iFluor™ 750 under optimal conditions
Application	Flow Cytometry (FACS)

**Spectral Properties**

---

Conjugate	PE/iFluor™ 750
Excitation Wavelength	566 nm
Emission Wavelength	778 nm

**Applications**

---

HIT11 is an anti-human monoclonal antibody that recognizes the CD2 antigen. CD2 (sometimes referred to as LFA-2) is a 45 kD single-pass type I membrane protein that is found on the surface of cells like T cells. CD2 has been thought to be involved with key biological processes such as cell-cell adhesion, especially heterotypic cell-cell adhesion. Also, in some organisms, it enhances myeloid dendritic cell activation, acts to positively regulate interferon-gamma secretion and is an enhancer of tumor necrosis factor production. CD2 is involved with key cellular pathways, for instance, the cell surface receptor signaling pathway. From a research standpoint, it is of biological interest due to its association

with critical macromolecules/ligands such as LFA-3, CD59, CD58 and CD48. CD2 is a moderately popular antibody target, with over 16000 publications in the last decade. CD2 is often used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of immunology. This antibody was purified through affinity chromatography and conjugated to PE/iFluor™ 750 (ex/em = 566/778 nm). It is compatible with the 561 nm laser and 783/56 nm bandpass filter (for example, as in the BD FACSMelody™).