

iFluor™ 405 Anti-human CD2 Antibody
HIT11Catalog number: 10020020, 10020021
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	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1
Immunogen	CD2 (LFA-2, Erythrocyte R, T11)
Clone	HIT11
Conjugate	iFluor™ 405

Biological Properties

Appearance	Light yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 405 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 405
Excitation Wavelength	403 nm
Emission Wavelength	427 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 iFluor™ 405 (ex/em = 403/427 nm). It is compatible with the 405 nm laser and 450/50 nm bandpass filter (for example, as in the BD FACSJazz™).