

**FITC Anti-rabbit/ dog/ mouse/ human
CD146 Antibody *P1H12***Catalog number: 114601G0, 114601G1
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	Rabbit, dog, mouse, human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1
Immunogen	CD146 (Muc-18, MCAM, Mel-CAM, S-endo)
Clone	P1H12
Conjugate	FITC

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with FITC under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	FITC
Excitation Wavelength	491 nm
Emission Wavelength	516 nm

Applications

P1H12 is an anti-rabbit/ dog/ mouse/ human monoclonal antibody that recognizes the CD146 antigen. CD146 (also known as MCAM) is a 118 kD member of the Ig superfamily that is expressed on the surface of cells such as T cells and endothelial cells. In some organisms, CD146 is involved in the positive regulation of cell migration, and is associated with a variety of biologically interesting macromolecules/ligands. CD146 is a fairly uncommon antibody target, with a little more than 4300 publications in the last decade. Even still, CD146 is often used in flow cytometry applications as a phenotypic marker for differentiation of cell types, specifically in the study of stem cells and innate immunity. This antibody

was purified through affinity chromatography and conjugated to FITC (ex/em = 491/516 nm). It is compatible with the 488 nm laser and 530/30 nm bandpass filter (for example, as in the BD FACSAria™ II).