

**iFluor™ 800 Anti-rabbit/ dog/ mouse/
human CD146 Antibody *P1H12***Catalog number: 114600N0, 114600N1
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	iFluor™ 800

Biological Properties

Appearance	Green liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 800 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 800
Excitation Wavelength	801 nm
Emission Wavelength	820 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 iFluor™ 800 (ex/em = 801/820 nm).