

**iFluor™ 860 Anti-mouse CD49 Antibody
*5H10-27 (MFR5)***Catalog number: 104930R0, 104930R1
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	Mouse
Class	Primary
Clonality	Monoclonal
Host	Rat
Isotype	Rat IgG2a kappa
Immunogen	CD49e (VLA-5 α , Integrin α 5 chain, ITGA5)
Clone	5H10-27 (MFR5)
Conjugate	iFluor™ 860

Biological Properties

Appearance	Green liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 860 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 860
Excitation Wavelength	853 nm
Emission Wavelength	878 nm

Applications

5H10-27 (MFR5) is an anti-mouse monoclonal antibody that forms an immune complex with the CD49e antigen. CD49e (also known as ITGA5, Integrin α 5 chain or VLA-5 α) is a 135 kD member of the Integrin family that is found on the surface of cells like stem cells and T cells. CD49 is associated with a variety of biologically interesting macromolecules/ligands, for example, fibronectin. CD49 is a relatively rare antibody target,

with fewer than 600 publications in the last decade. Even still, CD49e is commonly used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of cell biology. This antibody was purified through affinity chromatography and conjugated to iFluor™ 860 (ex/em = 853/878 nm).