

iFluor™ 405 Anti-human CD328 Antibody
6-434Catalog number: 13280020, 13280021
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	CD328 (Siglec7, AIRM1)
Clone	6-434
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

6-434 is an anti-human monoclonal antibody that forms an immune complex with the CD328 antigen. CD328 (alternatively called Siglec7 or AIRM1) is a 75 kD transmembrane protein that is found on the surface of cells like macrophages and T cells. CD328 is associated with a variety of biologically interesting macromolecules/ligands, in particular, silylated glycans. CD328 is a relatively rare antibody target, with fewer than 20

publications in the last decade. Even still, CD328 has a variety of applications in immunology research, often serving as a phenotypic marker for differentiating cell types in flow cytometric applications. 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 Miltenyi Biotec MACSQuant X).