

Purified Anti-human CD169 Antibody
7-239Catalog number: 11690000
Unit size: 100 ug**Product Details**

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	0.5 mg/mL
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide

Antibody Properties

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1
Immunogen	CD169 (Sialoadhesin, Siglec-1)
Clone	7-239

Biological Properties

Appearance	Liquid
Preparation	Antibody purified by affinity chromatography
Application	Flow Cytometry (FACS), ELISA, HC, Western Blot

Applications

7-239 is an anti-human monoclonal antibody that is specific for the CD169 antigen. CD169 (also known as Sialoadhesin) is a glycoprotein that is expressed on the surface of cells such as dendritic cells and macrophages. CD169 is associated with a variety of biologically interesting macromolecules/ligands, namely, 3-sialylated ligands, $\alpha 2$ and CD227. CD169 is a fairly uncommon antibody target, with a little more than 1900 publications in the last decade. Even still, CD169 is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of cell biology and immunology. This antibody was purified through affinity chromatography.