

iFluor™ 450 Anti-human CD169 Antibody
7-239Catalog number: 11690040, 11690041
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 | CD169 (Sialoadhesin, Siglec-1) |
| Clone | 7-239 |
| Conjugate | iFluor™ 450 |

Biological Properties

| | |
|-------------|--|
| Appearance | Brown liquid |
| Preparation | Antibody purified by affinity chromatography and then conjugated with iFluor™ 450 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging |

Spectral Properties

| | |
|-----------------------|-------------|
| Conjugate | iFluor™ 450 |
| Excitation Wavelength | 451 nm |
| Emission Wavelength | 502 nm |

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 and conjugated to iFluor™ 450 (ex/em = 451/502 nm). It is compatible with the 445 nm laser and 510/80 nm bandpass filter (for example, as in the BD FACSAria™ III).