

**mFluor™ Blue 570 Anti-human CD1
Antibody *HI149***Catalog number: 100100T0, 100100T1
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	CD1a (R4, T6)
Clone	HI149
Conjugate	mFluor™ Blue 570

Biological Properties

Appearance	Red liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ Blue 570 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	mFluor™ Blue 570
Excitation Wavelength	553 nm
Emission Wavelength	565 nm

Applications

HI149 is an anti-human monoclonal antibody that forms an immune complex with the CD1a antigen. CD1a (sometimes called R4 or T6) is a 49 kD member of the Ig superfamily that is found on the surface of cells like macrophages, dendritic cells and T cells. In certain organisms, CD1 is involved in the positive regulation of T cell mediated cytotoxicity, and is associated with a variety of biologically interesting

macromolecules/ligands, namely, β -2-Microglobulin. CD1 is a moderately popular antibody target, with over 15000 publications in the last decade. CD1a is typically used in flow cytometry applications as a phenotypic marker for differentiation of cell types, specifically in the study of immunology. This antibody was purified through affinity chromatography and conjugated to mFluor™ Blue 570 (ex/em = 553/565 nm). It is compatible with the 561 nm laser and 577/35 nm bandpass filter (for example, as in the Luminex Amnis ImageStream).