

PE/Texas Red® Anti-human CD1 Antibody
SN13Catalog number: 100121R0, 100121R1, 100121R2
Unit size: 25 tests, 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 kappa
Immunogen	CD1b (R1)
Clone	SN13
Conjugate	PE/Texas Red®

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/Texas Red® under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/Texas Red®
Excitation Wavelength	567 nm
Emission Wavelength	615 nm

Applications

The SN13 monoclonal antibody binds to human CD1b, a transmembrane protein commonly found on the surface of langerhans cells, dendritic cells and Tregs. In some organisms, CD1 is a positive regulator of T cell mediated cytotoxicity, and is associated with a variety of biologically interesting macromolecules/ligands, in particular, β -2-microglobulin. CD1 is a moderately popular antibody target, with over 15000 publications in the last decade. CD1b is commonly used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of immunology. This antibody was purified through affinity chromatography and conjugated to PE/Texas Red® (ex/em = 567/615 nm).

It is compatible with the 561 nm laser and 615/20 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte Advanteon).