

iFluor™ 710 Anti-human CD53 Antibody
HI29Catalog number: 105300K0, 105300K1
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	CD53 (Tetraspanin-25, MOX44)
Clone	HI29
Conjugate	iFluor™ 710

Biological Properties

Appearance	Blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 710 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 710
Excitation Wavelength	717 nm
Emission Wavelength	739 nm

Applications

HI29 is an anti-human monoclonal antibody that targets the CD53 antigen. CD53 (sometimes referred to as Tetraspanin-25, MOX44 or OX44) is a 35 - 42 kD member of the tetraspan family that is found on the surface of cells such as NK cells. In some organisms, CD53 enhances myoblast fusion, and is associated with a variety of biologically interesting macromolecules/ligands, in particular, VLA-4, Integrins and HLA-DR. CD53 is a

relatively rare antibody target, with fewer than 600 publications in the last decade. Even still, CD53 has a variety of applications in costimulatory molecules 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™ 710 (ex/em = 717/739 nm).