

**iFluor™ 514 Anti-human CD73 Antibody**  
**\*AD2\***Catalog number: 10730060, 10730061  
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 kappa
Immunogen	CD73 (NT5E, Ecto-5'-nucleotidase)
Clone	AD2
Conjugate	iFluor™ 514

**Biological Properties**

Appearance	Orange-red liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 514 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

**Spectral Properties**

Conjugate	iFluor™ 514
Excitation Wavelength	528 nm
Emission Wavelength	555 nm

**Applications**

AD2 is an anti-human monoclonal antibody that is specific for the CD73 antigen. CD73 (sometimes called NT5E or Ecto-5'-nucleotidase) is a 70 kD transmembrane protein that is found on the surface of cells such as epithelial cells, endothelial cells and stem cells. CD73 is associated with a variety of biologically interesting macromolecules/ligands, for instance, AMP. CD73 is a moderately popular antibody target, with over 11000

publications in the last decade. CD73 is commonly used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of stem cells. This antibody was purified through affinity chromatography and conjugated to iFluor™ 514 (ex/em = 528/555 nm). It is compatible with the 532 nm laser and 575/25 nm bandpass filter (for example, as in the Luminex Guava easyCyte).