

Purified Anti-human CD73 Antibody *AD2*

Catalog number: 10730000

Unit size: 100 ug

Product Details

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	0.5 mg/mL
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide

Antibody Properties

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1 kappa
Immunogen	CD73 (NT5E, Ecto-5'-nucleotidase)
Clone	AD2

Biological Properties

Appearance	Liquid
Preparation	Antibody purified by affinity chromatography
Application	Flow Cytometry (FACS), ELISA, HC, Western Blot

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.