

**PE/XFD610 Anti-human CD73 Antibody**  
**\*AD2, XFD610 Same Structure to Alexa**  
**Fluor™ 610\***Catalog number: 107301P0, 107301P1, 107301P2  
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	CD73 (NT5E, Ecto-5'-nucleotidase)
Clone	AD2
Conjugate	PE/AF610

**Biological Properties**

---

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/AF610 under optimal conditions
Application	Flow Cytometry (FACS)

**Spectral Properties**

---

Conjugate	PE/AF610
Excitation Wavelength	567 nm
Emission Wavelength	627 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 PE/XFD610 (ex/em =

567/627 nm). XFD610 is manufactured by AAT Bioquest, and it has the same chemical structure of Alexa Fluor® 610 (Alexa Fluor® is the trademark of ThermoFisher). It is compatible with the 561 nm laser and 615/20 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte).