

**iFluor™ 660 Anti-human CD39 Antibody**  
**\*TU66\***Catalog number: 103900G0, 103900G1  
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 IgG2b kappa
Immunogen	CD39 (ENTPD1, NTPDase 1)
Clone	TU66
Conjugate	iFluor™ 660

**Biological Properties**

Appearance	Blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 660 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

**Spectral Properties**

Conjugate	iFluor™ 660
Excitation Wavelength	663 nm
Emission Wavelength	678 nm

**Applications**

TU66 is an anti-human monoclonal antibody that recognizes the CD39 antigen. CD39 (also known as NTPDase 1 or ENTDP1) is a transmembrane protein that is found on the surface of cells such as dendritic cells and macrophages. CD39 is associated with a variety of biologically interesting macromolecules/ligands, for instance, ADP/ATP. CD39 is a fairly uncommon antibody target, with a little more than 4500 publications in the last

decade. Even still, CD39 is often 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 iFluor™ 660 (ex/em = 663/678 nm). It is compatible with the 640 nm laser and 670/30 nm bandpass filter (for example, as in the BD LSRFortessa™ X-20).