

mFluor™ UV460 Anti-human CD39 Antibody
TU66Catalog number: 103900Y0, 103900Y1
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	mFluor™ UV460

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ UV460 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	mFluor™ UV460
Excitation Wavelength	358 nm
Emission Wavelength	456 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 mFluor™ UV460 (ex/em = 358/456 nm). It

is compatible with the 355 nm laser and 450/50 nm bandpass filter (for example, as in the BD LSRFortessa™ Cell Analyzer).