

## XFD700 Anti-human CD30 Antibody \*MEM-268\*

Catalog Number: 103001A0, 103001A1

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	Lot specific (please consult certificate of analysis for given lot)
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 IgG
Immunogen	CD30 (Ber-H2, Ki-1, TNFRSF8)
Clone	MEM-268
Conjugate	AF700

### Biological Properties

---

Appearance	Dark blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF700 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging
Recommended Dilutions	For flow cytometry applications, the suggested concentration is at 5 uL/million cells in 100 uL staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined. <i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i>

### Spectral Properties

---

Conjugate AF700  
Excitation Wavelength 696 nm  
Emission Wavelength 719 nm

## Applications

---

MEM-268 is an anti-human monoclonal antibody that recognizes the CD30 antigen. CD30 (also known as Ber-H2, TNFRSF8 or Ki-1) is a 64 kD transmembrane protein that is located on the surface of cells like T cells and macrophages. In many organisms, CD30 promotes apoptotic process, is a suppressor of cell population proliferation and is an enhancer of TRAIL biosynthetic process. Additionally, it is a component of important cellular pathways, for instance, the tumor necrosis factor-mediated signaling pathway. From a research standpoint, it is of biological interest due to its association with essential macromolecules/ligands such as CD153. CD30 is a fairly uncommon antibody target, with a little more than 6900 publications in the last decade. Even still, CD30 is commonly used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of stem cells and immunology. This antibody was purified through affinity chromatography and conjugated to XFD700 (ex/em = 696/719 nm). XFD700 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 700 (Alexa Fluor® is the trademark of Thermo Fisher).