

Purified Mouse Anti-human/non-human primates Granzyme B Antibody *CLB-GB11, monoclonal*

Catalog number: V103760

Unit size: 0.1 mg

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	Tris buffered saline (TBS) solution, 15 mM sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human, non-human primates
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Granzyme B
Clone	CLB-GB11

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	FC (QC TESTED)

Applications

Granzyme B (alternatively called Cathepsin G-like 1, SECT, Human lymphocyte protein, Fragmentin-2, C11, HLP, CTSG1, T-cell serine protease 1-3E, Lymphocyte protease, Granzyme-2, CTLA-1 or Cytotoxic T-lymphocyte proteinase 2) is a protein with a molecular weight of 28 kDa, found in the nucleus, cytoplasm and cytosol of cells. In Homo sapiens, granzyme B is the subject of comprehensive research stemming from the fact that it plays a role in the positive regulation of protein insertion into mitochondrial membrane involved in apoptotic signaling pathway and granzyme-mediated apoptotic signaling pathway. Sequencing of granzyme B has shown it contains a primary structural unit, the peptidase S1 domain. Granzyme B has been thought to be involved with essential functions such as serine-type peptidase and serine-type endopeptidase activity. It aids in organismal processes, in particular, apoptotic process, natural killer cell mediated cytotoxicity and cytolysis. It acts to negatively regulate translation.