

**PE Mouse Anti-human/cow Perforin  
Antibody \*dG9, monoclonal, Cross  
Adsorbed\***

Catalog number: V1032005

Unit size: 100 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), 15 mM sodium azide, 0.2% (w/v) BSA

**Antibody Properties**

---

Species Reactivity	Human, cow
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Perforin
Clone	dG9
Conjugate	PE

**Biological Properties**

---

Preparation	Antibody purified by affinity chromatography, cross-adsorbed against mouse serum and then conjugated with PE under optimal conditions
Application	FC (QC TESTED)

**Applications**

---

Perforin-1 (sometimes called Cytolysin, Lymphocyte pore-forming protein or PFP) is a protein with a molecular weight of 61 kDa, expressed in the membrane, cytolytic granule and extracellular region of cells. In humans, perforin-1 is involved in the positive regulation of killing of cells of other organism. Perforin-1 recognizes calcium ion and identical protein, and additionally, takes part in organismal processes, in particular, immunological synapse formation, protein homooligomerization and apoptotic process. Sequencing of perforin-1 has demonstrated it contains 3 types of conserved structural units: EGF-like, C2 and MACPF domain. Perforin-1 has been closely linked to vital functions such as wide pore channel activity. It is clinically significant because abnormalities in its function have been associated with diseases like familial hemophagocytic lymphohistiocytosis 2 (FHL2). Familial hemophagocytic lymphohistiocytosis 2, an autosomal recessive inheritance disorder characterized by hepatomegaly, hypofibrinogenemia and lymphadenopathy, has in particular been of interest to researchers.