

**Purified Mouse Anti-human HLA-E Antibody**  
**\*MEM-E/07, monoclonal\***

Catalog number: V1031060

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	Phosphate-buffered saline (PBS, pH 7.2), 15 mM sodium azide, 0.2% (w/v) BSA

**Antibody Properties**

---

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	HLA-E
Clone	MEM-E/07

**Biological Properties**

---

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

**Applications**

---

HLA class I histocompatibility antigen,  $\alpha$  chain E is a 40 kDa transmembrane protein that can be found in the plasma membrane, cell surface and extracellular space of cells. Sequencing of HLA class I histocompatibility antigen,  $\alpha$  chain E has shown it contains 3 types of conserved structural units: extracellular, cytoplasmic and Ig-like C1-type domain. HLA class I histocompatibility antigen,  $\alpha$  chain E takes part in processes such as immune response and natural killer cell mediated immunity. It promotes antibody-dependent cellular cytotoxicity, tumor necrosis factor production and TRAIL production conversely also inhibits natural killer cell mediated cytotoxicity. HLA class I histocompatibility antigen,  $\alpha$  chain E is the subject of comprehensive study in part because of the fact that it is involved with the type I interferon signaling pathway and interferon- $\gamma$ -mediated signaling pathway. HLA class I histocompatibility antigen,  $\alpha$  chain E binds to MHC class I protein, signaling receptor and T cell receptor.