

# Purified Low Endotoxin Mouse Anti-human/cow/non-human primates HLA Class I Antibody \*W6/32, monoclonal, Cross Adsorbed\*

Catalog number: V103850

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	Low endotoxin azide free Phosphate-buffered saline (PBS, pH 7.2), 0.2 µm filter sterilized. Endotoxin level is less than 0.01 EU/µg of the protein, as determined by the LAL test., 0.2% (w/v) BSA

## Antibody Properties

---

Species Reactivity	Human, cow, non-human primates, cat
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	HLA Class I
Clone	W6/32
Conjugate	Purified Low Endotoxin

## Biological Properties

---

Preparation	Antibody purified by affinity chromatography, cross-adsorbed against rabbit serum and then conjugated with Purified Low Endotoxin under optimal conditions
Application	FC (QC TESTED), IP, WB, IHC(F), ICC, ELISA, MC (CYTOF), FUNC

## Applications

---

HLA class I histocompatibility antigen,  $\alpha$  chain G is a 44 kDa transmembrane protein that can be expressed in the plasma membrane, integral component of lumenal side of endoplasmic reticulum membrane and MHC class I protein complex of cells. It is also known as MHC class I antigen G and HLA G antigen. In Homo sapiens, HLA class I histocompatibility antigen,  $\alpha$  chain G plays an important role in organismal processes, in particular, cellular defense response, antigen processing and presentation of endogenous peptide antigen via MHC class Ib and protection from natural killer cell mediated cytotoxicity. It recognizes peptide antigen, signaling receptor and CD8 receptor. HLA class I histocompatibility antigen,  $\alpha$  chain G is an enhancer of T cell tolerance induction, cellular senescence and macrophage cytokine production in contrast to also is an inhibitor of protein kinase B signaling, G0 to G1 transition and immune response. It has been thought to be involved with critical functions like protein homodimerization activity. Sequencing of HLA class I histocompatibility antigen,  $\alpha$  chain G has exemplified it contains 3 conserved structural units: extracellular, cytoplasmic and Ig-like C1-type domain. HLA class I histocompatibility antigen,  $\alpha$  chain G is the subject of extensive study stemming from the fact that it plays a role in the type I interferon signaling pathway, immune response-inhibiting cell surface

receptor signaling pathway and interferon- $\gamma$ -mediated signaling pathway. HLA class I histocompatibility antigen,  $\alpha$  chain G is an integral part of immune response.