

**Purified Mouse Anti-human/mouse/rabbit  
MHC Class I ( $\alpha$ + $\beta$ 2m chains) Antibody**  
**\*IVA26, monoclonal\***

Catalog number: V1031755  
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, mouse, rabbit, pig, cow, sheep, coturnix japonica |
| Class              | Primary                                                  |
| Clonality          | Monoclonal                                               |
| Host               | Mouse                                                    |
| Immunogen          | MHC Class I ( $\alpha$ + $\beta$ 2m chains)              |
| Clone              | IVA26                                                    |

### Biological Properties

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

### Applications

Major histocompatibility complex class I-related gene protein is a transmembrane protein with a molecular weight of 45 kDa, found in the endoplasmic reticulum membrane, late endosome membrane and endoplasmic reticulum of cells. In humans, major histocompatibility complex class I-related gene protein has been closely linked to important functions like MHC class I receptor activity. It binds with  $\beta$ -2-microglobulin and T cell receptor. Sequencing of major histocompatibility complex class I-related gene protein has exemplified it contains 3 conserved structural units: extracellular, cytoplasmic and Ig-like C1-type domain. It acts to positively regulate T cell mediated cytotoxicity directed against tumor cell target. It aids in organismal processes, for instance, immune response, T cell differentiation in thymus and cytokine production involved in immune response.