

**Description**

The MAGE-A1 Peptide (278-286, KVLEYVIKV) is a peptide corresponding to MAGE-A1 (Melanoma-associated antigen 1), amino acids 278-286. The MAGE-A1 peptide is widely used to stimulate human MAGE-A1-specific CD8<sup>+</sup> T cells or as negative control in the stimulation of human MAGE-A4-specific CD8<sup>+</sup> T cells.

**Background**

MAGE (melanoma associated antigen) proteins are CT (cancer testis) antigens, and there are about 60 proteins in the MAGE family that can be subdivided into type I (present only on the X-chromosome, MAGE-A, B, and C) and type II (MAGE D-L and necdin). Under normal conditions they are mostly found in the testis and placenta. They are found at high levels in several cancer types, such as melanoma, brain, and breast cancer, and are involved in the development of resistance to chemotherapy, cell motility and cell survival. Expression of MAGE proteins tend to correlate with a poor prognosis. They are intracellular proteins, with MAGE-A1 being mostly cytosolic, making them poor targets for strategies such as CAR-T cell therapy. MAGE proteins are degraded in the proteasome, and the peptides created can then be found on the cell membrane in combination with MHC (major histocompatibility complex) I. The presentation on the cell surface in this form makes them an attractive target for TCR (T cell receptor)-T cell therapy. Several clinical trials are ongoing, and either alone or in combination with other forms of cancer therapy the use of TCR-T cells targeting MAGE-A antigens may become a promising therapeutical avenue.

**Sequence**

KVLEYVIKV

**Species**

Human

**Supplied As**

Liquid, 100 µl

**Formulation**

1 mM peptide in DMSO

**Stability**

At least one year at -80°C.

**Storage**

Upon first thaw, aliquot and store at -80°C. Avoid repeated freeze-thaw cycles.

**Application**

Stimulation of human MAGE-A1-specific CD8<sup>+</sup> T cells or as negative control in the stimulation of human MAGE-A4-specific CD8<sup>+</sup> T cells.

**Related Products**

<i>Products</i>	<i>Catalog #</i>	<i>Size</i>
MAGE-A1-Specific TCR Lentivirus	78934	2 x 500 µl
MAGE-A4 Peptide (230-239)	78966	100 µl

Version 022824