

Description

The Melanoma-associated antigen recognized by T cells-1 (MART-1) peptide (26-35, Leu27; ELAGIGILTV) is a peptide corresponding to MART-1 amino acids 26-35 in which the Alanine (A) was replaced with Leucine (L) at position 27. This peptide has higher binding affinity towards MART-1-specific T cells and is more stable than the native MART-1 peptide (26-35, EAAGIGILTV; BPS Bioscience #78759).

Background

MART-1 (also known as Melan-A) is a differentiation antigen expressed at the surface of melanocytes. A peptide fragment of the protein is found bound to MHC complexes that are recognized by cytotoxic T cells. MART-1 is found in most skin cancers, including melanomas, and is used as a biomarker for diagnostic purposes. Since the expression of the protein is restricted to melanocytes containing tissues (skin and retina) and is not found in other tissues, MART-1 is an attractive target for cancer vaccines and adoptive cell therapy. MART-1 peptide 26-35 is a fragment commonly associated with MHC and recognized by T cell receptors.

Sequence

ELAGIGILTV

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 MART-1-specific CD8⁺ T cells.

Related Products

<i>Products</i>	<i>Catalog #</i>	<i>Size</i>
NY-ESO-1-Specific TCR Lentivirus (Clone c259)	78676	100 µl/2 x 500 µl
NY-ESO-1-Specific TCR Lentivirus (Clone 1G4)	78675	100 µl/2 x 500 µl
MART-1-Specific TCR Lentivirus (Clone DMF4)	78678	100 µl/2 x 500 µl
MART-1-Specific TCR Lentivirus (Clone DMF5)	78679	100 µl/2 x 500 µl
MART-1 peptide (27-35)	78761	100 µl
MART-1 peptide (26-35)	78759	100 µl
NY-ESO-1 Peptide (157-165)	78758	100 µl