

Instruction Manual

Catalog Number	PK-AB577-5823
Description	MART-1 (also known as Melan-A) is a melanocyte differentiation antigen recognized by autologous cytotoxic T lymphocytes. Six other melanoma associated antigens recognized by autologous cytotoxic T cells include MAGE-1, Tyrosinase, gp100, gp75, BAGE-1, and GAGE-1. Subcellular fractionation shows that MART-1 is present in melanosomes and endoplasmic reticulum.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	E. coli expressed recombinant human MART-1
Clone / IgG Subtype	Rabbit IgG
Species Reactivity	human
Specificity	See Applications or Species Reactivity.
Formulation	100 µg (0.5 mg/ml) affinity purified rabbit polyclonal antibody in phosphate-buffered saline (PBS) containing 30% glycerol, 0.5% BSA, and 0.01% thimerosal.
Reconstitution	Can be diluted in other aqueous buffers at the concentrations determined for the respective application.
Storage & Stability	Store at -20°C. For long-term storage, aliquot and refreeze at -70°C. Avoid repeated freeze/thaw cycles.
Applications	Western blotting (0.5-4 µg/ml). The optimal concentrations should be determined individually. Antibody might be suitable for other applications not tested so far.

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.