

MART-1 Antibody [#C0257]

- Catalog Number:** C0257-1, C0257-2, C0257
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** Q16655
- All Names:** MART-1, MART1, melan-A, MELAN A, MELANOMA ANTIGEN RECOGNIZED BY T CELLS 1
- Form of Antibody:** Rabbit IgG in phosphate buffered saline (without Mg²⁺ and Ca²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
- Storage/Stability:** Store at -20°C/1 year
- Immunogen:** The antiserum was produced against synthesized peptide derived from human MART-1.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** MART-1 antibody detects endogenous levels of total MART-1 protein.
- Reactivity:** Human
- Applications:** WB: 1:300~1:1000 IHC: 1:50~1:100 ELISA: 1:40000
- References:** Toshihiko Hoashi, J. Biol. Chem., Apr 2005; 280: 14006 - 14016.
Maria P. Bettinotti, J. Immunol., Jul 1998; 161: 877.
P. J. Frost, J. Immunol., Mar 2001; 166: 3564.
Flora Wang, Clin. Cancer Res., Oct 1999; 5: 2756.