

## MCL1 Antibody [#C0258]

- Catalog Number:** C0258-1, C0258-2, C0258
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** Q07820
- All Names:** EAT/MCL-1 protein, Induced myeloid leukemia cell differentiation protein Mcl-1, Myeloid cell leukemia sequence 1.
- Form of Antibody:** Rabbit IgG in phosphate buffered saline (without Mg<sup>2+</sup> and Ca<sup>2+</sup>), 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 MCL1.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** MCL1 antibody detects endogenous levels of total MCL1 protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000
- References:** Alfredo De Biasio, J. Biol. Chem., Jun 2007; 10.1074.  
Potchanapond Graidist, J. Biol. Chem., Sep 2004; 279: 40868 - 40875.  
Kenichi Fujise, J. Biol. Chem., Dec 2000; 275: 39458.  
Di Zhang, J. Biol. Chem., Sep 2002; 277: 37430 - 37438.