

HDAC1 Antibody [#C0221]

- Catalog Number:** C0221-1, C0221-2, C0221
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No.:** Q13547
- All Names:** Histone deacetylase 1, HD1, RPD3L1
- 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 HDAC1.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** HDAC1 antibody detects endogenous levels of total HDAC1 protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:300~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
- References:** Richard L. Cai, J. Biol. Chem., Sep 2000; 275: 27909.
Takayoshi Matsumura, J. Biol. Chem., Apr 2005; 280: 12123 - 12129.
Gregory David, J. Biol. Chem., Jun 2002; 277: 23658 - 23663.
Yuh-Long Chang, PNAS, Aug 2001; 98: 9730.