

HDAC2 (Ab-394) Antibody [#B7099]

- Catalog Number:** B7099-1, B7099-2, B7099
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- 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 non-phosphopeptide derived from human HDAC2 around the phosphorylation site of Serine 394.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity:** HDAC2(Ab-394) Antibody detects endogenous levels of total HDAC2 protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100
ELISA: 1:1000
- Swiss-Prot No. :** Q92769
- References:** Tsai SC, et al.(2002)J Biol Chem; 277(35): 31826-33.