

## JAB1 Antibody [#C0237]

- Catalog Number:** C0237-1, C0237-2, C0237
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No.:** Q9UNS2
- All Names:** SGN3, COP9, SUBUNIT 3
- 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 JAB1.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** JAB1 antibody detects endogenous levels of total JAB1 protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:5000
- References:** George Z. Rassidakis, Clin. Cancer Res., Mar 2003; 9: 1121.  
Ricardo V. Lloyd, Am. J. Pathol., Mar 2002; 160: 1171.  
Shu Xiao, PNAS, Oct 2006; 103: 16448 - 16453.  
Alex Kudrin, J. Biol. Chem., Oct 2006; 281: 29641 - 29651.