

## Ret (Ab-905) Antibody [#B7213]

- Catalog Number:** B7213-1, B7213-2, B7213
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** P07949
- All Names:** Amplexin, CTTN, EMS1, Oncogene EMS1, SRC9, Src substrate cortactin
- 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 non-phosphopeptide derived from human Ret around the phosphorylation site of tyrosine 905 (K-L-Y<sup>P</sup>-G-M).
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity:** Ret (Ab-905) antibody detects endogenous levels of total Cortactin protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1:1000 ELISA: 1:10000
- References:** Domenico Salvatore, Cancer Res., Feb 2001; 61: 1426 - 1431.  
Hye Sook Jung, Mol. Endocrinol., Nov 2005; 19: 2748 - 2759.  
Iván Plaza-Menacho, J. Biol. Chem., Mar 2007; 282: 6415 - 6424.  
Matthias Drosten, J Natl Cancer Inst, Aug 2004; 96: 1231 - 1239.