



## Caveolin-1(Ab-14) Antibody

Catalog Number 21112

(For Research Use Only)

- Catalog Number:** 21112-1, 21112-2, 21112  
**Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl  
**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 Caveolin-1 around the phosphorylation site of Tyrosine 14.  
**Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.  
**Specificity/Sensitivity:** Caveolin-1(Ab-14) Antibody detects endogenous levels of total Caveolin-1 protein.  
**Reactivity:** Human, Mouse, Rat  
**Applications:** WB: 1:500~1:1000 ELISA: 1:20000  
**Swiss-Prot No. :** Q03135  
**References:** Zhang Y, et al. (2005) Mol Cell Proteomics. 4(9): 1240-1250.  
Labrecque L, et al. (2004) J Biol Chem. 279(50): 52132-52140.  
Fielding PE, et al. (2004) Biochemistry. 43(9): 2578-2586.  
Labrecque L, et al. (2003) Mol Biol Cell. 14(1): 334-347.  
Maggi D, et al. (2002) Biochem Biophys Res Commun. 295(5): 1085-1089.

