



## Connexin43 (Ab-367) Antibody

Catalog Number 21250

(For Research Use Only)

- Catalog Number:** 21250-1, 21250-2, 21250  
**Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl  
**Swiss-Prot No. :** P17302  
**All Names:** Cx43; Gap junction 43 kDa heart protein  
**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 phosphopeptide derived from human Connexin43 around the phosphorylation site of serine 367 (R-A-S<sup>P</sup>-S-R).  
**Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.  
**Specificity/Sensitivity** Connexin43 (Ab-367) antibody detects endogenous levels of total Connexin43 protein.  
**Reactivity:** Human, Mouse, Rat  
**Applications:** WB: 1:500~1:1000 ELISA: 1:20000  
**References:** Xiaoyong Bao, et al. (2004) Am J Physiol Cell Physiol; 286: C647-C654.  
K Boengler, et al. (2006) Heart; 92: 1724-1727.  
Satoshi Matsushita, et al. (2006) J. Histochem. Cytochem; 54: 343-353.  
Holger K. Eltzschig, et al. (2006) Circ. Res; 99: 1100-1108.

