



Connexin43 (Phospho-Ser367) Antibody

Catalog Number 11258

(For Research Use Only)

- Catalog Number:** 11258-1, 11258-2, 11258
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** P17302
- All Names:** Cx43; Gap junction 43 kDa heart protein
- All Sites:** Human: Ser367; Mouse: Ser367; Rat: Ser367
- 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 phosphopeptide derived from human Connexin43 around the phosphorylation site of serine 367 (R-A-S^P-S-R).
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific phosphopeptide. The antibody against non-phosphopeptide was removed by chromatography using non-phosphopeptide corresponding to the phosphorylation site.
- Specificity/Sensitivity:** Connexin43 (Phospho-Ser367) antibody detects endogenous levels of Connexin43 only when phosphorylated at serine 367.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100
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.

