

Cox1 Antibody [#C0159]

- Catalog Number:** C0159-1, C0159-2, C0159
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** P23219
- All Names:** COX3, PCOX1, CYCLOOXYGENASE 1, PARTIAL COX1 PROTEINS, PGHS1
CYCLOOXYGENASE 3, PROSTAGLANDIN G/H SYNTHASE 1, prostaglandin-
endoperoxide synthase 1
- 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 peptide derived from human COX1.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using
epitope-specific immunogen.
- Specificity/Sensitivity** COX1 antibody detects endogenous levels of total COX1 protein.
- Reactivity:** Human
- Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:40000
- References:** Lawrence M. Szewczuk, J. Biol. Chem., May 2004; 279: 22727 - 22737.
Syed R. Baber, Am J Physiol Heart Circ Physiol, Oct 2005; 289: H1476 - H1487.
Subhashini Chandrasekharan, J. Lipid Res., Dec 2005; 46: 2636 - 2648.
K. Trent Moreland, J. Pharmacol. Exp. Ther., May 2007; 321: 446 - 454.