

PTGR1 Antibody [#C16862]

Catalog Number:	C16862
Concentration:	1mg/ml
Swiss-Prot No.:	Q14914
Other Names:	Prostaglandin reductase 1; PRG-1; NADP-dependent leukotriene B4 12-hydroxydehydrogenase; EC=1.3.1.74; 15-oxoprostaglandin 13-reductase; EC=1.3.1.48; LTB4DH
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg ²⁺ and Ca ²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized peptide derived from internal of human PTGR1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PTGR1 antibody detects endogenous levels of total PTGR1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Humphray S.J., Nature 429:369-374(2004). The MGC Project Team; Genome Res. 14:2121-2127(2004). Yokomizo T., J. Biol. Chem. 271:2844-2850(1996).