

SCN4B Antibody [#C18808]

Catalog Number:	C18808
Concentration:	1mg/ml
Swiss-Prot No.:	P16949
Other Names:	Sodium channel subunit beta-4 [Precursor]; SCN4B
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 SCN4B.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	SCN4B antibody detects endogenous levels of total SCN4B protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Yu F.H., J. Neurosci. 23:7577-7585(2003). The MGC Project Team; Genome Res. 14:2121-2127(2004). Medeiros-Domingo A., Circulation 116:134-142(2007).