

## OMG Antibody [#C17547]

<b>Catalog Number:</b>	C17547
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P23515
<b>Other Names:</b>	Oligodendrocyte-myelin glycoprotein
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	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.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human OMG.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	OMG antibody detects endogenous levels of total OMG protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100 ELISA: 1:10000
<b>References:</b>	Viskochil D., Mol. Cell. Biol. 11:906-912(1991). Mikol D.D., J. Cell Biol. 110:471-479(1990). Mikol D.D., J. Cell Biol. 111:2673-2679(1990).