

## hnRNP G Antibody [#C10581]

<b>Catalog Number:</b>	C10581
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P38159
<b>Other Names:</b>	HNRPG, RBMX, RNA binding motif protein, X chromosome, ROG, glycoprotein p43, heterogeneous nuclear ribonucleoprotein G
<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 N-terminal of human hnRNP G.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	hnRNP G antibody detects endogenous levels of total hnRNP G protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:5000
<b>References:</b>	Soulard M., Nucleic Acids Res. 21:4210-4217(1993). Ross M.T., Nature 434:325-337(2005). Venables J.P., Hum. Mol. Genet. 8:959-969(1999).