

## PMP22 Antibody [#C0306]

<b>Catalog Number:</b>	C0306
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q01453
<b>Other Names:</b>	GAS3, GROWTH ARREST-SPECIFIC 3
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without $Mg^{2+}$ and $Ca^{2+}$ ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from C-terminal of human PMP22.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	PMP22 antibody detects endogenous levels of total PMP22 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:40000
<b>References:</b>	GJ Snipes, J. Cell Biol., Apr 1992; 117: 225. P A Csurhes, J. Neurol. Neurosurg. Psychiatry, Oct 2005; 76: 1431 - 1439. Laura D. Attardi, Genes & Dev., Mar 2000; 14: 704. Julie A. Hodapp, Arch Neurol, Jan 2006; 63: 112 - 117.