

## AGR3 Antibody [#C11390]

<b>Catalog Number:</b>	C11390
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q8TD06
<b>Other Names:</b>	AG-3; hAG-3; AG3; Breast cancer membrane protein 11
<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 AGR3.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	AGR3 antibody detects endogenous levels of total AGR3 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
<b>References:</b>	Adam P.J., J. Biol. Chem. 278:6482-6489(2003). Clark H.F., Genome Res. 13:2265-2270(2003). Zhang Z., Protein Sci. 13:2819-2824(2004).