

## C3AR1 Antibody [#G028]

<b>Catalog Number:</b>	G028
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q16581
<b>Other Names:</b>	C3a anaphylatoxin chemotactic receptor; C3a-R; C3AR; C3AR1; AZ3B, C3R1, HNFAG09
<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 C3AR1.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	C3AR1 antibody detects endogenous levels of total C3AR1 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
<b>References:</b>	Roglic A., Biochim. Biophys. Acta 1305:39-43(1996). Crass T., Eur. J. Immunol. 26:1944-1950(1996). Ames R.S., J. Biol. Chem. 271:20231-20234(1996).