

## ADORA3 Antibody [#G204]

<b>Catalog Number:</b>	G204
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P33765
<b>Other Names:</b>	Adenosine A3 receptor;ADORA3; UNQ1931/PRO4406
<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 ADORA3.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	ADORA3 antibody detects endogenous levels of total ADORA3 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 92%, Positives = 92%); Rat (Identities = 92%, Positives = 92%)
<b>Applications:</b>	IF: 1:500~1:1000 ELISA: 1:10000
<b>References:</b>	Sajjadi F.G., <i>Biochim. Biophys. Acta</i> 1179:105-107(1993). Salvatore C.A., <i>Proc. Natl. Acad. Sci. U.S.A.</i> 90:10365-10369(1993). Meyerehof W., Submitted (DEC-1993) to the EMBL/GenBank/DDBJ databases.