

## OLFML2A Antibody [#G484]

<b>Catalog Number:</b>	G484
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q68BL7
<b>Other Names:</b>	olfactomedin-like; olfactomedin-like 2A; OLFML2A; photomedin-1
<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 OLFML2A.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	OLFML2A antibody detects endogenous levels of total OLFML2A protein.
<b>Reactivity:</b>	Human
<b>Applications:</b>	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:10000
<b>References:</b>	Clark H.F., Genome Res. 13:2265-2270(2003). Humphray S.J., Nature 429:369-374(2004). The MGC Project Team; Genome Res. 14:2121-2127(2004).