

## STAM2 (Ab-192) Antibody [#B1237]

<b>Catalog Number:</b>	B1237
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	O75886
<b>Other Names:</b>	HBP, HRS binding protein, Signal transducing adaptor molecule (SH3 domain and ITAM motif) 2, Signal transducing adaptor molecule 2A, Signal transducing adaptor molecule 2B
<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 non-phosphopeptide derived from human STAM2 around the phosphorylation site of tyrosine 192 (S-L-Y <sup>P</sup> -P-S).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	STAM2 (Ab-192) antibody detects endogenous levels of total STAM2 protein.
<b>Reactivity:</b>	Human, Mouse
<b>Applications:</b>	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:5000
<b>References:</b>	Endo K., FEBS Lett. 477:55-61(2000). Wiemann S., Genome Res. 11:422-435(2001). Hillier L.W., Nature 434:724-731(2005).