

## Dab1 (Ab-220) Antibody [#B7054]

<b>Catalog Number:</b>	B7054
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	O75553
<b>Other Names:</b>	disabled 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 non-phosphopeptide derived from human Dab1 around the phosphorylation site of tyrosine 220 (N-I-Y <sup>P</sup> -Q-V).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	Dab1 (Ab-220) antibody detects endogenous levels of total Dab1 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:100 IF: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Lambert de Rouvroit C., Genomics 53:246-247(1998). Fazili Z., Submitted (MAY-2000) to the EMBL/GenBank/DDBJ databases. Gregory S.G., Nature 441:315-321(2006).