

## CD71/TfR (Ab-24) Antibody [#B0852]

<b>Catalog Number:</b>	B0852
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P02786
<b>Other Names:</b>	CD71 antigen, T9, TFRC, TR, TfR1, Transferrin receptor protein 1, Trfr, p90
<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 CD71/TfR around the phosphorylation site of serine 24 (R-F-S <sup>P</sup> -L-A).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	CD71/TfR (Ab-24) antibody detects endogenous levels of total CD71/TfR protein.
<b>Reactivity:</b>	Human, Mouse
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:10000
<b>References:</b>	Evans P., Gene 199:123-131(1997). Gene 199:123-131(1997). Coppolino M., Biochem. J. 306:129-134(1995). Liu T., J. Proteome Res. 4:2070-2080(2005).