

AhR (Ab-36) Antibody [#B0765]

Catalog Number:	B0765
Concentration:	1mg/ml
Swiss-Prot No.:	P35869
Other Names:	Ah receptor, Aryl hydrocarbon receptor precursor
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg^{2+} and Ca^{2+}), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized non-phosphopeptide derived from human AhR around the phosphorylation site of serine 36 (N-P-S ^P -K-R).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	AhR (Ab-36) antibody detects endogenous levels of total AhR protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:100~1:500 ELISA: 1:5000
References:	Itoh S., Nucleic Acids Res. 21:3578-3578(1993). Dolwick K.M., Mol. Pharmacol. 44:911-917(1993). Hillier L.W., Nature 424:157-164(2003).