

Androgen Receptor (Ab-650) Antibody [#B7009]

Catalog Number:	B7009
Concentration:	1mg/ml
Swiss-Prot No.:	P10275
Other Names:	ANDR; Androgen receptor; DHTR; Dihydrotestosterone receptor; NR3C4
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 Androgen Receptor around the phosphorylation site of serine 650 (T-T-S ^P -P-T).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Androgen Receptor (Ab-650) antibody detects endogenous levels of total Androgen Receptor protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:40000
References:	Lubahn D.B., Mol. Endocrinol. 2:1265-1275(1988). Lubahn D.B., Proc. Natl. Acad. Sci. U.S.A. 86:9534-9538(1989). Govindan M.V.; Mol. Endocrinol. 4:417-427(1990).