Androgen Receptor (Ab-363) Antibody [#B8204]

Catalog Number:B8204Concentration:1mg/mlSwiss-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 Mg²⁺ and Ca²⁺), 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 tyrosine 363 (D-Y-Y^P-N-F).

Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-

specific immunogen.

Specificity: Androgen Receptor (Ab-363) antibody detects endogenous levels of total Androgen Receptor

protein.

Reactivity: Human (Identities = 100%, Positives = 100%); Mouse (Identities = 92%, Positives = 100%);

Rat (Identities = 92%, Positives = 100%)

Applications: WB: 1:500~1:1000 ELISA: 1:10000

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).

Technical Support: support@aatbio.com; 408-733-1055