

AR

Androgen Receptor

human, recombinant, Sf9 Insect cells

Cat. No.	Amount
PR-867	5 µg

For *in vitro* use only
Quality guaranteed for 12 months
Store at -80°C

Avoid freeze / thaw cycles

Form

Liquid. Supplied in 20 mM Tris-Cl, 25% Glycerol, 100 mM KCl, 1 mM DTT, 0.2 mM EDTA

Application

Use only for research and not for drug or diagnostic purposes.

AR can be used for gel mobility shift assay, for protein-protein and small molecules-protein interactions assay.

Purity

> 95% by SDS-PAGE.

Description

AR (dihydrotestosterone receptor; testicular feminization; spinal and bulbar muscular atrophy; Kennedy disease). The androgen receptor is an androgen-activated member of the nuclear receptor superfamily of transcription factors. Signaling by the androgen receptor plays a key role in proper development and function of male reproductive organs. The androgen receptor gene is more than 90 kb long and codes for a protein that has 3 major functional domains: the N-terminal domain, DNA-binding domain, and androgen-binding domain. The protein functions as a steroid-hormone activated transcription factor. Upon binding the hormone ligand, the receptor dissociates from accessory proteins, translocates into the nucleus, dimerizes, and then stimulates transcription of androgen responsive genes. This gene contains 2 polymorphic trinucleotide repeat segments that encode polyglutamine and polyglycine tracts in the N-terminal transactivation domain of its protein. Expansion of the polyglutamine tract causes spinal bulbar muscular atrophy (Kennedy disease). Mutations in this gene are also associated with complete androgen insensitivity (CAIS). Two alternatively spliced variants encoding distinct isoforms have been described. The androgen receptor (AR) has been shown to play a critical role in the development and progression of the prostate cancer.

Selected References:

- Marques, R.B. et al. (2005) Androgen receptor modifications in prostate cancer cells upon long-term androgen ablation and antiandrogen treatment. *Int. J. Cancer* **117**(2):221-229.
- Metzger, E. et al. (2005) LSD1 demethylates repressive histone marks to promote androgen-receptor-dependent transcription. *Nature* **437**(7057):436-439.
- Murthy, S. et al. (2005) Androgen receptor signaling and vitamin D receptor action in prostate cancer cells. *Prostate* **64**(4):362-372.
- Hillmer, A.M. et al. (2005) Genetic variation in the human androgen receptor gene is the major determinant of common early-onset androgenetic alopecia. *Am. J. Hum. Genet.* **77**(1):140-148.
- Gao, S. et al. (2005) The androgen receptor directly targets the cellular Fas/FasL-associated death domain protein-like inhibitory protein gene to promote the androgen-independent growth of prostate cancer cells. *Mol. Endocrinol.* **19**(7):1792-1802.
- Chen, G. et al. (2005) Androgen receptor mutants detected in recurrent prostate cancer exhibit diverse functional characteristics. *Prostate* **63**(4):395-406.
- Lim, J.T. et al. (2005) Cyclin-dependent kinase 6 associates with the androgen receptor and enhances its transcriptional activity in prostate cancer cells. *Proc. Natl. Acad. Sci. USA* **102**(14):5156-5161.
- Kinoshita, M. et al. (2005) Differently regulated androgen receptor transcriptional complex in prostate cancer compared with normal prostate. *Am. J. Hum. Genet.* **12**(4):390-397.
- Estebanez-Perpina, E. et al. (2005) The molecular mechanisms of

AR

Androgen Receptor

human, recombinant, Sf9 Insect cells

coactivator utilization in ligand-dependent transactivation by the androgen receptor. *J. Biol. Chem.* **280(9)**:8060-8068.