Androgen receptor monoclonal antibody (SPM335)

The expression of AR is reportedly inversely correlated with histologic grade i.e. well differentiated prostate tumors show higher expression than the poorly differentiated tumors. In prostate cancer, AR has been proposed, as a marker of hormone-responsiveness and thus it may be useful in identifying patients likely to benefit from anti-androgen therapy. Anti-androgen receptor has been useful clinically in differentiating morpheaform basal cell carcinoma (mBCC) from desmoplastic trichoepithelioma (DTE) in the skin.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

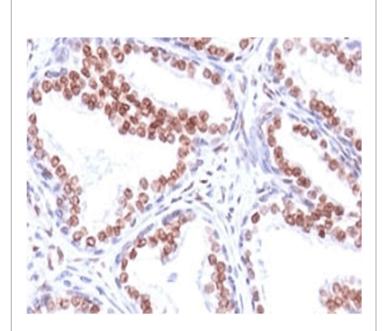
Order Online »

ENZ-ABS770-0100

100µg

Manuals, SDS & CofA

View Online »



IHC: Formalin-fixed, paraffin-embedded human prostate carcinoma stained with anti-Androgen Receptor antibody

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Dihydrotestosterone receptor, Nuclear receptor subfamily

3 group C member 4

Application IHC

Application NotesThis mAb is superb for staining of formalin/paraffin tissues.

Clone SPM335

Crossreactivity Does not cross react with estrogen, progesterone, or

glucocorticoid receptors.

Formulation Liquid. In 1X PBS containing 0.1 mg/ml BSA and 0.05%

sodium azide.

Host Mouse

Immunogen Human androgen receptor, amino acids 302-318

(STEDTAEYSPFKGGYTK).

Isotype IgG1ĸ

Purity Detail Protein G affinity purified.

Species Reactivity Human

Specificity Recognizes a protein of 110kDa, which is identified as

androgen receptor (AR). It reacts with full length, and the

newly described A form of the receptor.

UniProt ID P10275

Last modified: May 29, 2024



info-

eu@enzolifesciences.com