

# P504S + p63 + HMWCK antibody cocktail

## Ready-to-use antibody cocktail to P504S (AMACR), p63 and high molecular weight

**Cytokeratins for use with IHC applications.**  
P504S, also known as  $\alpha$ -methylacyl coenzyme A racemase (AMACR), is a peroxisomal and mitochondrial enzyme that plays a role in bile acid synthesis and  $\beta$ -oxidation of branched chain fatty acids. P504S, expression has been found in prostatic adenocarcinoma and prostate glands involved in prostatic intraepithelial neoplasia, but not in benign glands or normal prostatic tissue. p63, a member of the p53 family of transcription factors, has been identified in nuclei of the basal epithelium in normal prostate glands and prostatic intraepithelial neoplasia, but not seen expressed in the vast majority of adenocarcinomas of the prostate. High molecular weight cytokeratins (HMWCK) CK5 and CK14 are expressed in a variety of normal and neoplastic epithelial tissues such as basal epithelia in the prostate gland.

This antibody is covered by our [Worry-Free Guarantee](#).

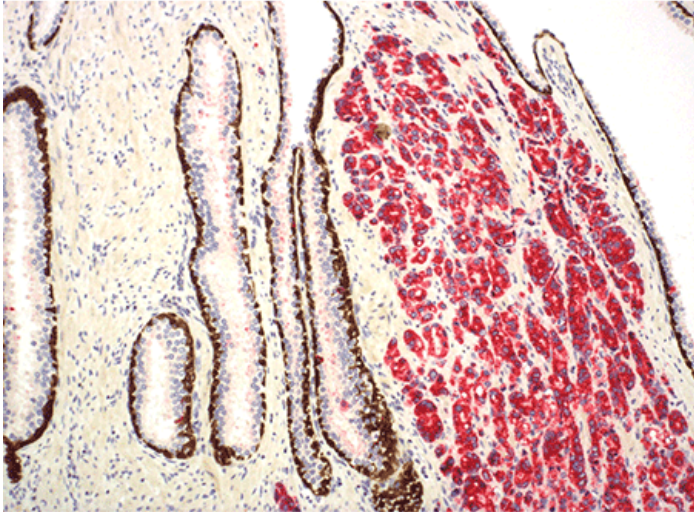
### Ordering Information

[Order Online »](#)

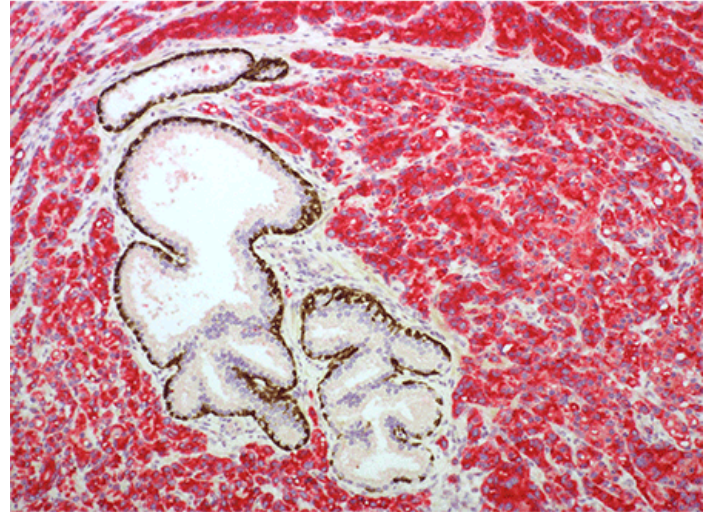
ENZ-ABS717-0010	10ml
-----------------	------

### Manuals, SDS & CofA

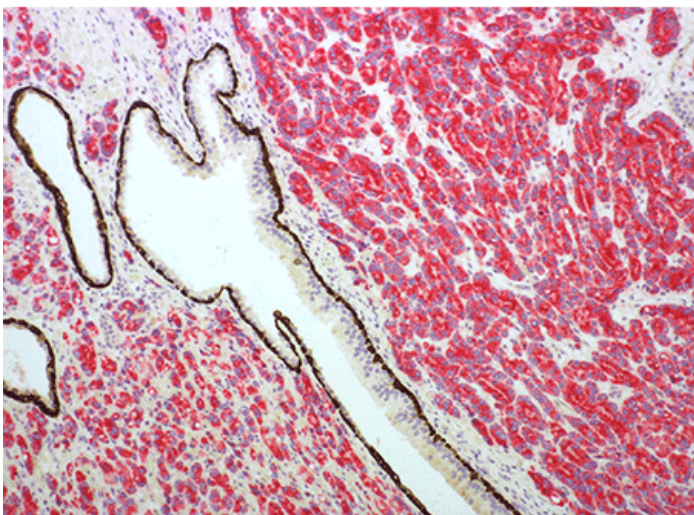
[View Online »](#)



Prostate tissue incubated with Enzo's P504S + p63 + HMWCK antibody cocktail (Prod. No. ENZ-ABS717). IHC staining was performed on an automated Leica BOND III. P504 stains (cytoplasm red) p63 (nuclear brown) and high molecular weight cytokeratins (cytoplasmic brown).



Formalin-fixed paraffin-embedded prostate tissue stained using Enzo's P504S+p63+HMWCK cocktail antibody (Prod. No. ENZ-ABS717). IHC staining was performed on an automated Leica BOND III. P504 stains (cytoplasm red) p63 (nuclear brown) and high molecular weight cytokeratins (cytoplasmic brown).



Formalin-fixed paraffin-embedded prostate tissue stained using Enzo's P504S+p63+HMWCK cocktail antibody (Prod. No. ENZ-ABS717). IHC staining was performed on an automated Leica BOND III. P504 stains (cytoplasm red) p63 (nuclear brown) and high molecular weight cytokeratins (cytoplasmic brown).

## Handling & Storage

Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	+4°C
Shipping	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Application	IHC
Application Notes	The P504S + p63 + HMWCK antibody cocktail can be used to qualitatively identify the presence of cytokeratin 1, 5, 10 & 14 (Cytoplasmic brown), p63 (nuclear brown) and P504S (cytoplasmic red) proteins in FFPE tissue sections using IHC test methods.
Formulation	Liquid. In DPBS buffer, pH 7.4, containing 1% BSA and 0.05% preservative.
Host	Rabbit, Mouse, Mouse
Immunogen	A peptide corresponding to residues in human P504S (AMACR), recombinant p63, keratin solubilized from human stratum corneum.
Isotype	IgG
Purity Detail	Affinity-purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (RTU)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	Q9UHK6 (P504S), Q9H3D4 (p63), P04264 (CK1), P13647 (CK5), P13645 (CK10), P02533 (CK14)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)