

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 α -methylacyl coenzyme A racemase (AMACR), is a peroxisomal and mitochondrial enzyme that plays a role in bile acid synthesis and β -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](#).

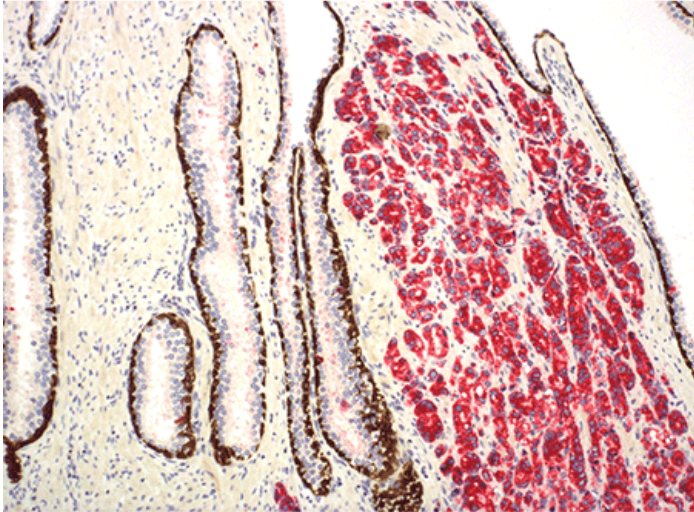
Ordering Information

[Order Online »](#)

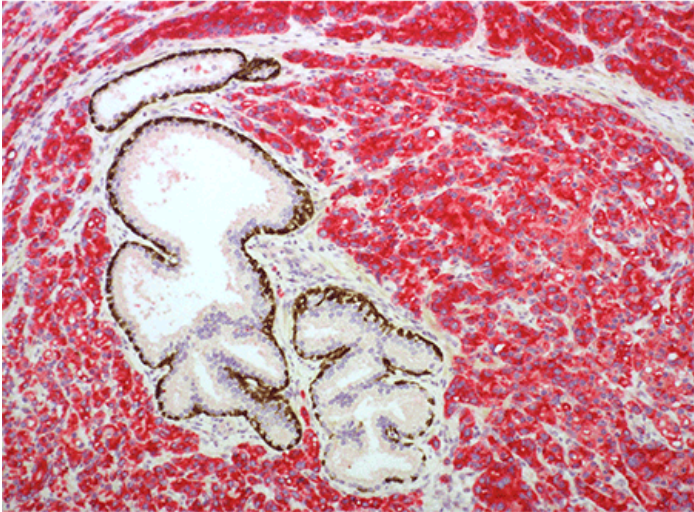
| | |
|-----------------|------|
| ENZ-ABS717-0010 | 10ml |
| ENZ-ABS717-0025 | 25ml |

Manuals, SDS & CofA

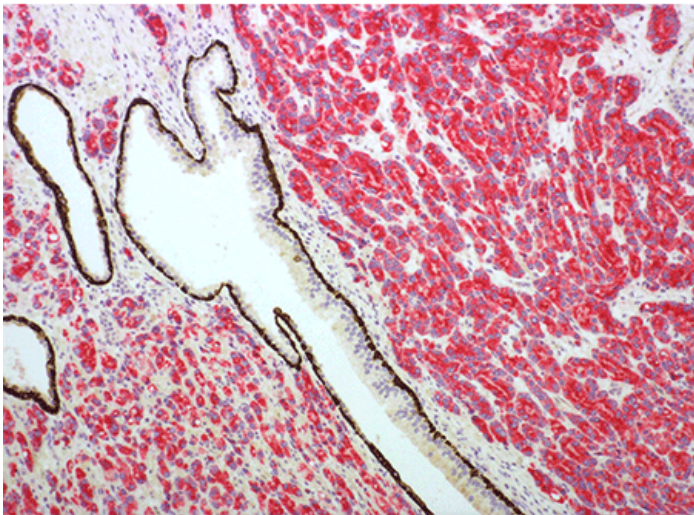
[View Online »](#)



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) |

Worry-free Guarantee

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

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)