

NKX3.1 monoclonal antibody (361)

NKX3.1 is a prostate-specific homeobox gene and transcription factor. It plays a key role in prostate development and tumor suppression, and is a reliable marker for prostate cancer in tissue analysis.

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

Ordering Information

[Order Online »](#)

ENZ-ABS1090-0500	500µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Formalin fixed paraffin embedded human prostate
stained with NKX3.1 antibody

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Homeobox protein Nkx-3.1

Application IHC (PS)

Application Notes Pretreatment – EDTA Buffer pH 8.0. Incubation & Temperature 30 min @ RT

Clone 361

Formulation Tissue culture supernatant. Contains sodium azide as a preservative.

Host Mouse

Immunogen Recombinant His fusion protein corresponding to full length human NKX3.1.

Isotype IgG1κ

Positive Control Prostate Cancer

Recommendation Dilutions/Conditions Immunohistochemistry (paraffin sections) (1:50-1:100)

Suggested dilutions/conditions may not be available for all applications.

Optimal conditions must be determined individually for each application.

Species Reactivity Human

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



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)