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

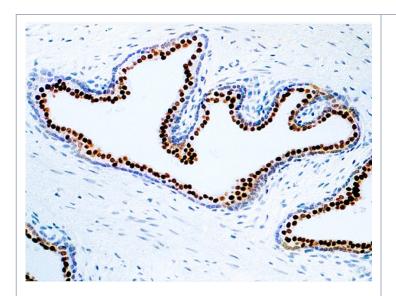
Order Online »

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.

lgG1κ

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

•

Last modified: June 6, 2025



info-

eu@enzolifesciences.com