

Kallikrein-3

monoclonal antibody

(03)

This mouse monoclonal antibody is specifically designed for the detection of KLK3 (Kallikrein-Related Peptidase 3), commonly known as prostate-specific antigen (PSA).

Kallikrein-3, also known as prostate-specific antigen (PSA), is a serine protease predominantly produced by the epithelial cells of the prostate gland. PSA plays a vital role in semen liquefaction by cleaving gel-forming proteins, facilitating sperm motility. Beyond its physiological function, PSA is a well-established biomarker for prostate cancer, benign prostatic hyperplasia (BPH), and other prostate-related conditions. Elevated PSA levels in the blood are commonly associated with prostate malignancies, making it a critical target for both diagnostic and therapeutic research in urology and oncology.

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

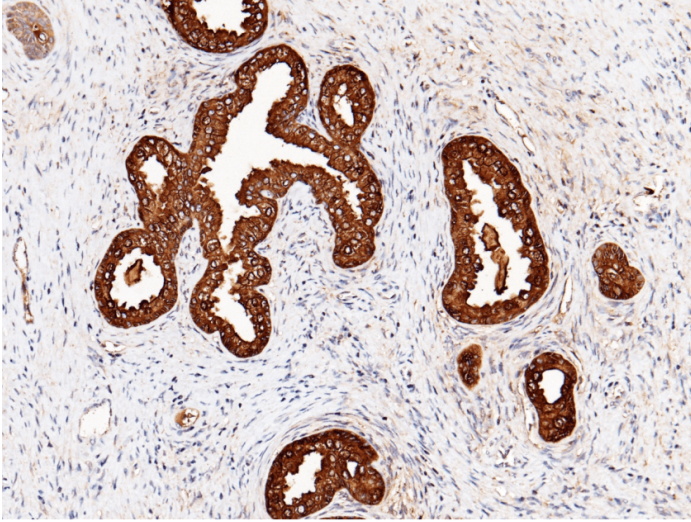
Ordering Information

[Order Online »](#)

ENZ-ABS1048-0100	100µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemical staining of human KLK3 in human prostate with mouse monoclonal antibody (1:60, formalin-fixed paraffin embedded sections).

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	PSA, Prostate-specific antigen
Application	IHC (PS)
Clone	03
Formulation	Liquid. 0.2 µm filtered solution in PBS. Carrier-free. Preservative-free.
Host	Mouse
Immunogen	Recombinant human Kallikrein-3 protein.
Isotype	IgG
Purity Detail	Protein A purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (paraffin sections) (1:50-1:200) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.
Source	This antibody was produced from a hybridoma resulting from the fusion of a mouse myeloma with B cells obtained from a mouse immunized with purified, recombinant human Kallikrein-3.
Species Reactivity	Human

Worry-free Guarantee

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

Last modified: January 9, 2025



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)