

Pan Cytokeratin polyclonal antibody

Cytokeratins are intermediate filament proteins found in epithelial cells. Pan cytokeratin antibodies recognize multiple cytokeratin subtypes, making them ideal for broadly identifying epithelial cell populations and epithelial-derived tumors in spatial studies.

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

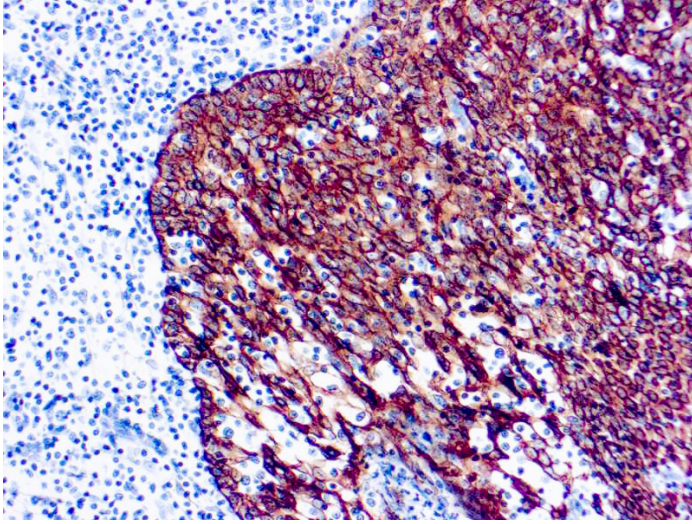
Ordering Information

[Order Online »](#)

ENZ-ABS1081-0500	500µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Formalin fixed paraffin embedded human colon adenocarcinoma stained with Pan Cytokeratin 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 pan-CK

Application IHC (PS)

Formulation Contains sodium azide as a preservative.

Host Rabbit

Immunogen Recombinant full-length human KRT76 and KRT77 proteins.

Isotype IgG

Positive Control Adeno Colon

Purity Detail Purified immunoglobulin.

Recommendation Immunohistochemistry (paraffin sections) (1:50-1:200)

Dilutions/Conditions 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)