

PAX8 monoclonal antibody

Monoclonal antibody to PAX-8 for use with IHC applications

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

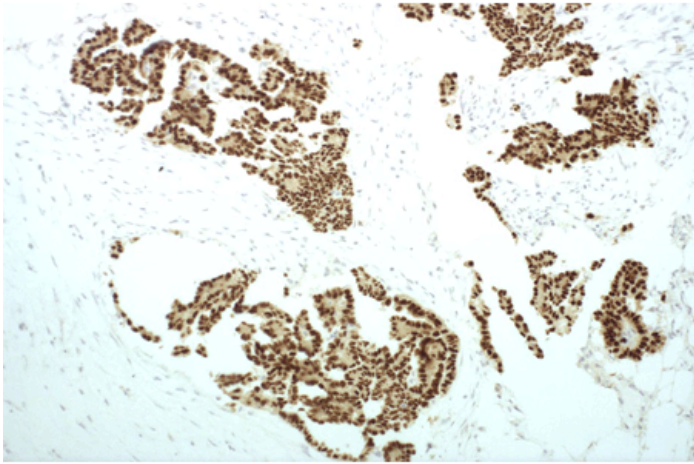
Ordering Information

[Order Online »](#)

ENZ-ABS718-1000	1ml
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Formalin-fixed paraffin-embedded serous ovarian carcinoma tissue stained using Enzo's PAX-8 antibody (Prod. No. ENZ-ABS718).

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

Alternative Name	Paired box protein Pax-8
Application	IHC
Application Notes	For IHC applications: Dilute the concentrate as appropriate in antibody diluent, and then apply 150 µL to the slide, as indicated in the user’s IHC protocol.
Formulation	Liquid. In Tris buffer, pH 7.5, containing 1% BSA and <0.1% sodium azide.
Host	Mouse
Immunogen	Recombinant PAX-8.
Isotype	IgG
Purity Detail	Affinity-purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (1:50-1:200)Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	Q06710
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)