

SARS2 polyclonal antibody

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

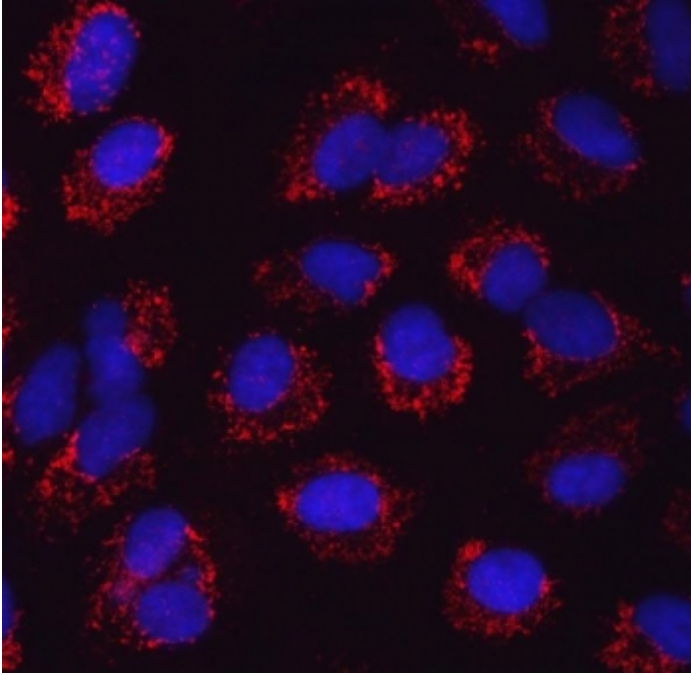
Ordering Information

[Order Online »](#)

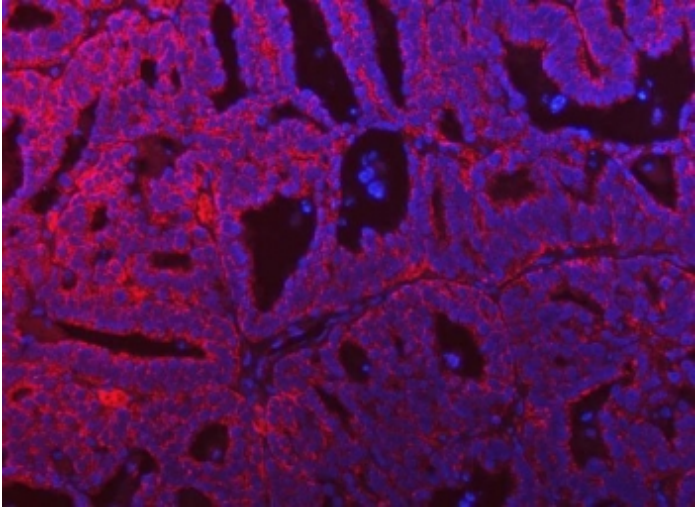
ENZ-ABS919-0100	100µg
-----------------	-------

Manuals, SDS & CofA

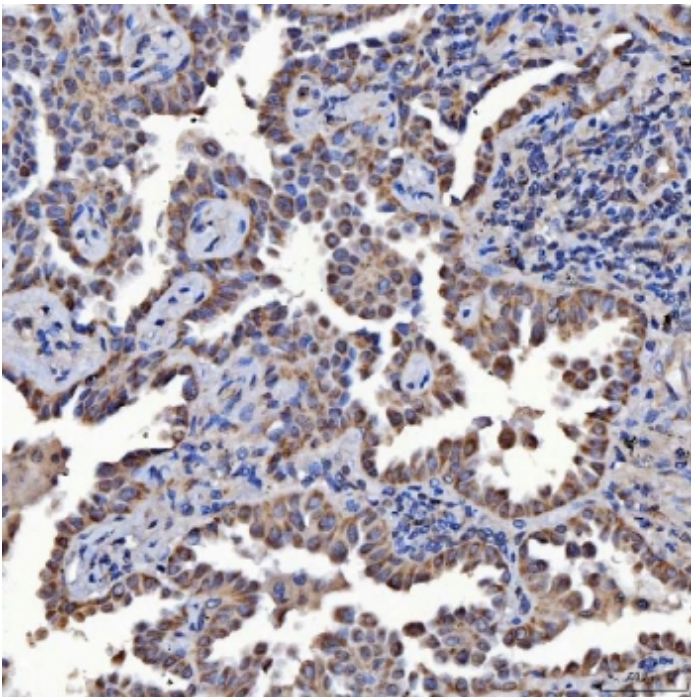
[View Online »](#)



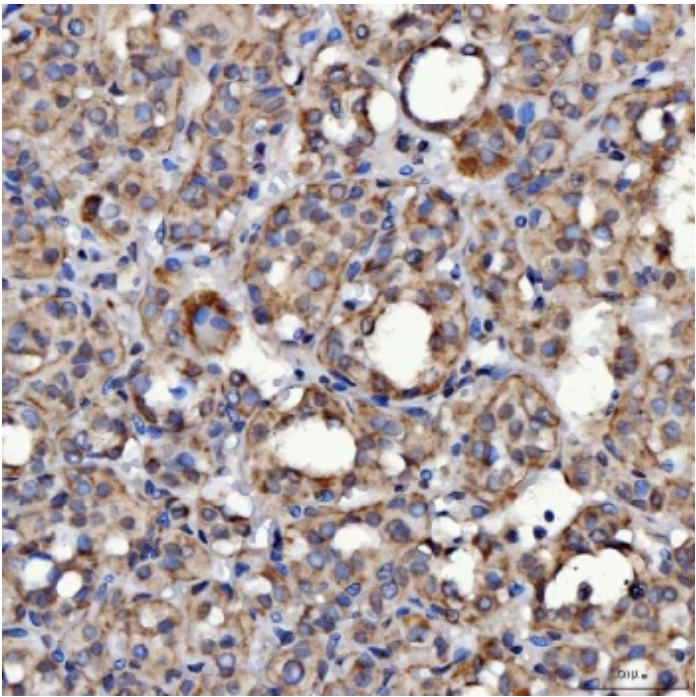
Immunofluorescent staining of FFPE human A549 cells with SerRS antibody (red) and DAPI nuclear stain (blue). HIER: steam section in pH6 citrate buffer for 20 min.



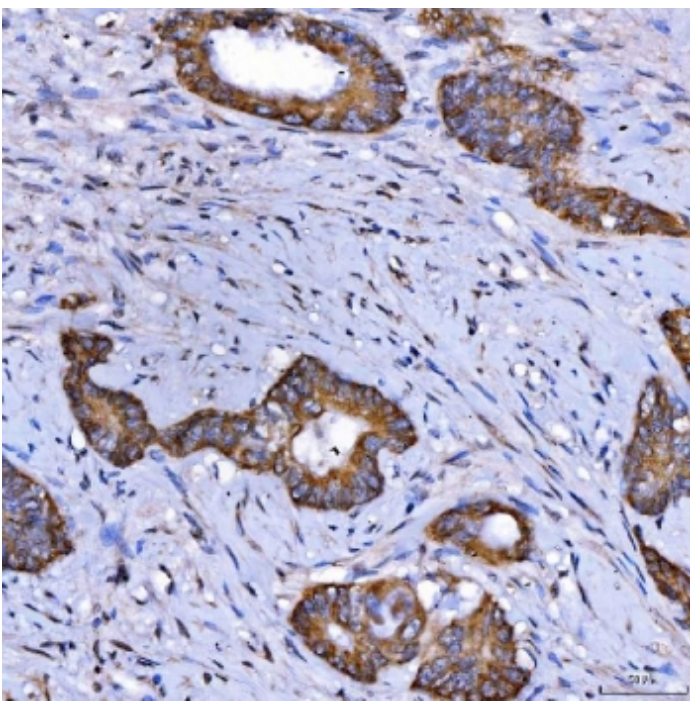
Immunofluorescent staining of FFPE human ovarian cancer tissue with SerRS antibody (red) and DAPI nuclear stain (blue). HIER: steam section in pH8 EDTA buffer for 20 min.



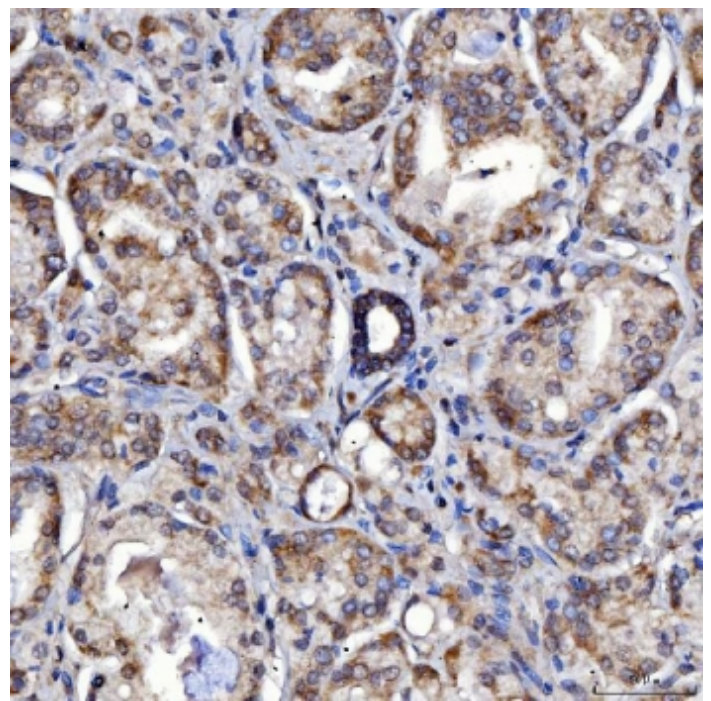
IHC staining of FFPE human lung cancer tissue with SerRS antibody. HIER: boil tissue sections in pH8 EDTA for 20 min and allow to cool before testing.



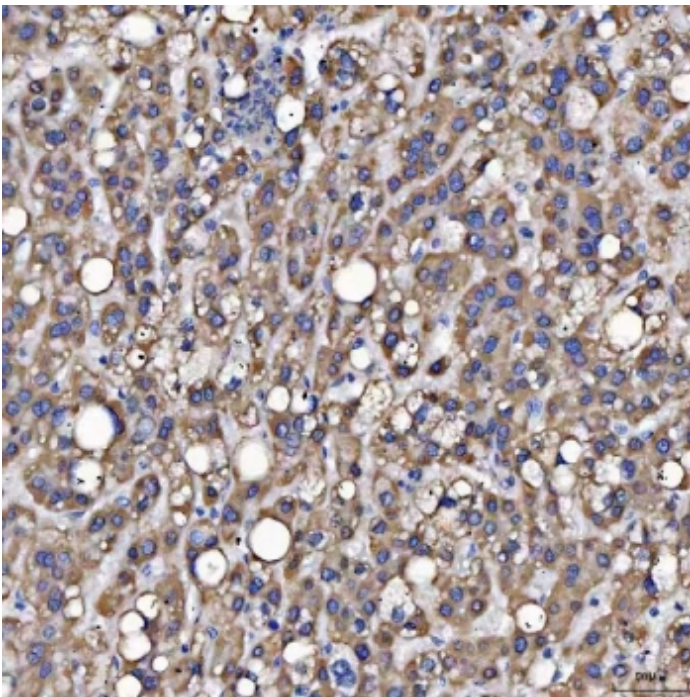
IHC staining of FFPE human thyroid cancer tissue with SerRS antibody. HIER: boil tissue sections in pH8 EDTA for 20 min and allow to cool before testing.



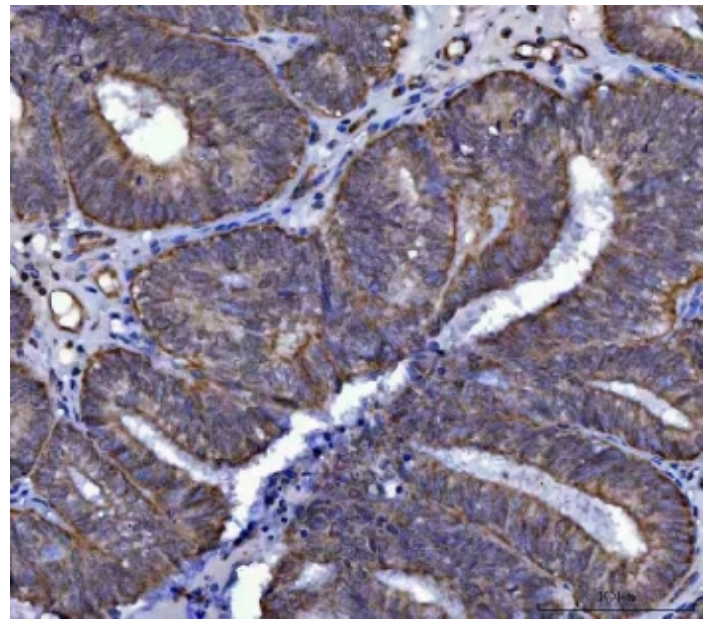
IHC staining of FFPE human colon cancer tissue with SerRS antibody. HIER: boil tissue sections in pH8 EDTA for 20 min and allow to cool before testing.



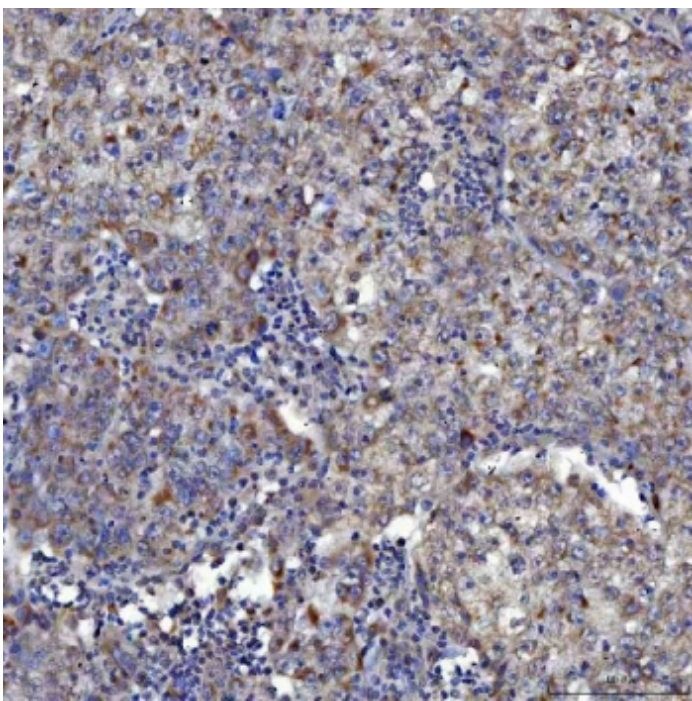
IHC staining of FFPE human prostate adenocarcinoma tissue with SerRS antibody. HIER: boil tissue sections in pH8 EDTA for 20 min and allow to cool before testing.



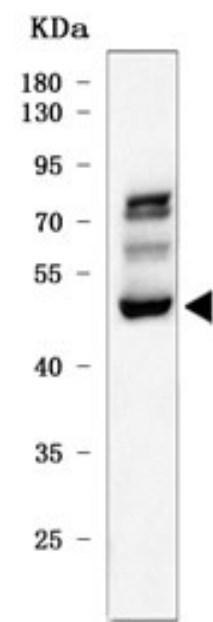
IHC staining of FFPE human liver cancer tissue with SerRS antibody. HIER: boil tissue sections in pH8 EDTA for 20 min and allow to cool before testing.



IHC staining of FFPE human endometrial cancer tissue with SerRS antibody. HIER: boil tissue sections in pH8 EDTA for 20 min and allow to cool before testing.



IHC staining of FFPE human cervical cancer tissue with SerRS antibody. HIER: boil tissue sections in pH8 EDTA for 20 min and allow to cool before testing.



Western blot testing of human HepG2 cell lysate with SerRS antibody. Predicted molecular weight ~58 kDa, commonly observed at 50-60 kDa.

Handling & Storage

Use/Stability	After reconstitution, the SerRS antibody can be stored for up to one month at +4°C. For long-term, aliquot and store at -20°C.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	SerRS, Seryl-tRNA synthetase
Application	ELISA, IF, IHC (PS), WB
Formulation	Lyophilized from 1X PBS containing 2% Trehalose
Host	Rabbit
Immunogen	E. coli-derived recombinant human SARS2 (aa 46-518).
Isotype	IgG
Purity Detail	Antigen affinity purified.
Recommendation Dilutions/Conditions	ELISA (direct) (0.1-0.5µg/ml)Immunofluorescence (5µg/ml)Immunohistochemistry (FFPE) (2-5µg/ml)Western blot (0.5-1µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	Q9NP81

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-
fr@enzolifesciences.com](mailto:info-fr@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)