

# S100A8 recombinant monoclonal antibody (216)

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

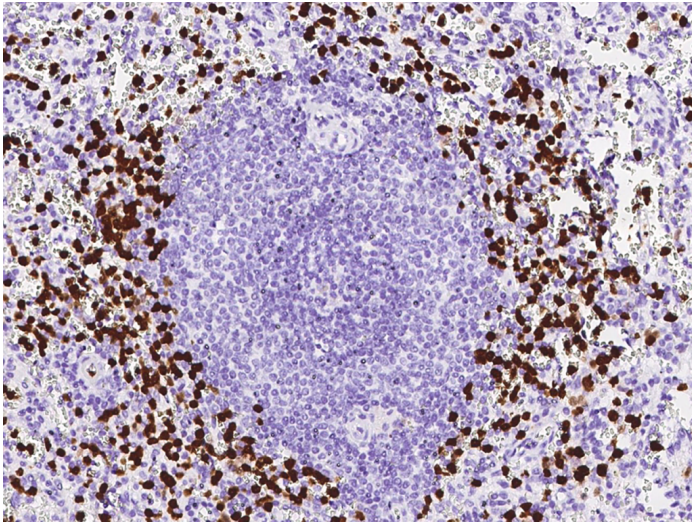
## Ordering Information

[Order Online »](#)

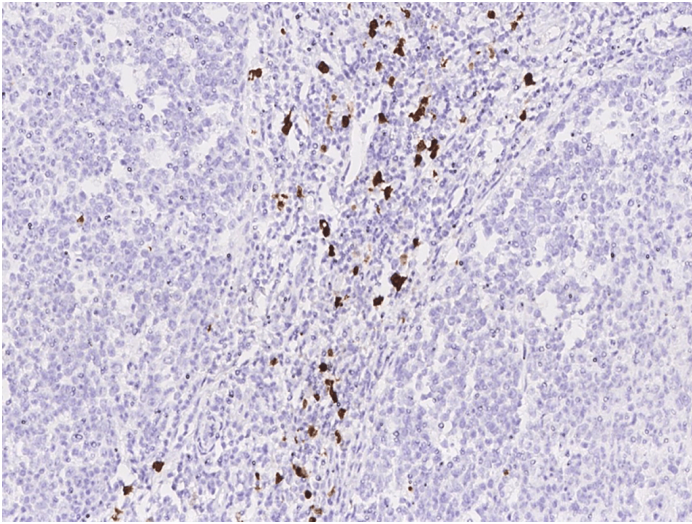
|                 |       |
|-----------------|-------|
| ENZ-ABS973-0100 | 100µl |
|-----------------|-------|

## Manuals, SDS & CofA

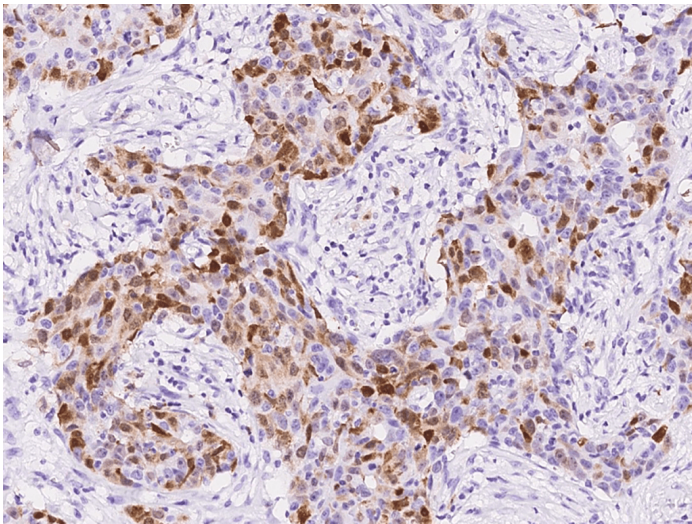
[View Online »](#)



Immunochemical staining of human S100A8 in human spleen with rabbit monoclonal antibody at 1:2000 dilution, formalin-fixed paraffin embedded sections.



Immunochemical staining of human S100A8 in human tonsil with rabbit monoclonal antibody at 1:2000 dilution, formalin-fixed paraffin embedded sections.



Immunochemical staining of human S100A8 in human breast carcinoma with rabbit monoclonal antibody at 1:2000 dilution, formalin-fixed paraffin embedded sections.

## Handling & Storage

**Handling** Protected from light. Avoid freeze-thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Application** IHC (PS)

**Clone** 216

**Formulation** Liquid. 0.2 µm filtered solution in PBS. Preservative free.

**Host** Rabbit

**Immunogen** Recombinant human S100A8 protein.

**Isotype** IgG

**Purity Detail** Protein A purified.

**Recommendation Dilutions/Conditions** Immunohistochemistry (paraffin sections) (1:500-1:5'000) 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: September 17, 2024**



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)