

# SLAMF7 monoclonal antibody (SLAMF7/3649)

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

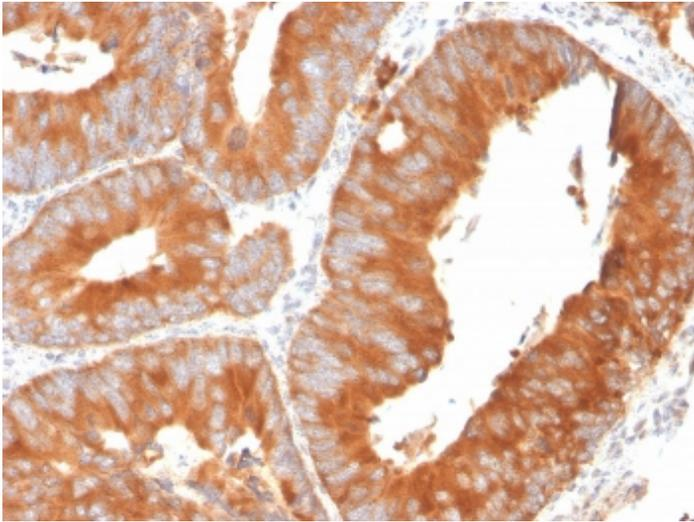
## Ordering Information

[Order Online »](#)

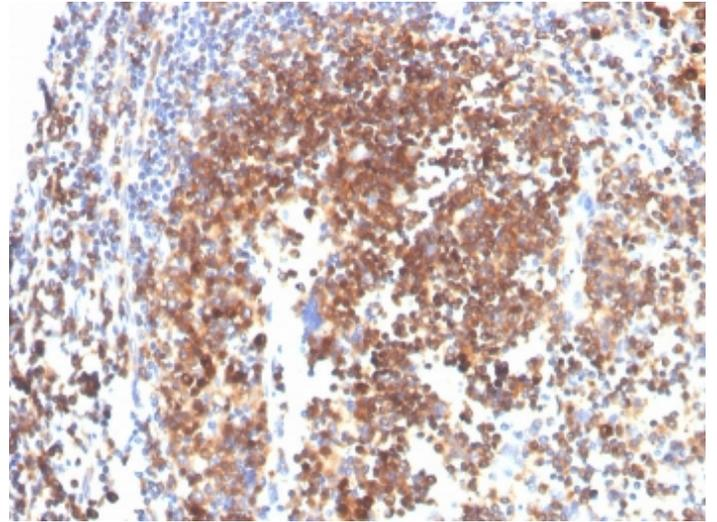
ENZ-ABS920-0100	100µg
-----------------	-------

## Manuals, SDS & CofA

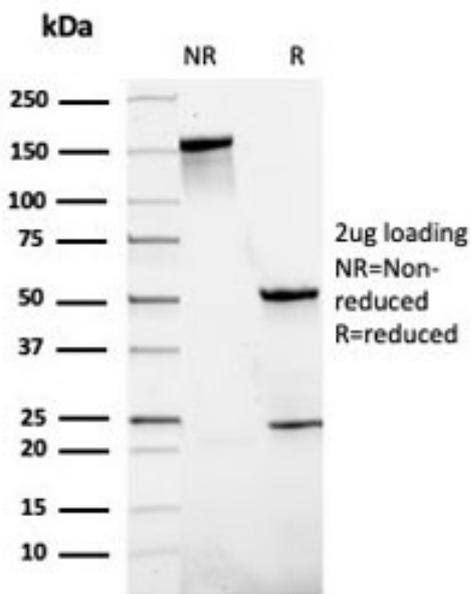
[View Online »](#)



IHC staining of FFPE human spleen with SLAMF7 antibody. HIER: boil tissue sections in pH 9 10mM Tris with 1mM EDTA for 20 min and allow to cool before testing.



IHC staining of FFPE human tonsil with SLAMF7 antibody. HIER: boil tissue sections in pH 9 10mM Tris with 1mM EDTA for 20 min and allow to cool before testing.



SDS-PAGE analysis of purified, BSA-free SLAMF7 antibody as confirmation of integrity and purity.

## Handling & Storage

<b>Handling</b>	Store the SLAMF7 antibody at 2-8°C (with azide) or aliquot and store at -20°C or colder (without azide).
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	CS1, CD319
<b>Application</b>	IHC (PS)
<b>Clone</b>	SLAMF7/3649
<b>Formulation</b>	Liquid. 0.2 mg/ml in 1X PBS containing 0.1 mg/ml BSA (US sourced), 0.05% sodium azide.
<b>Host</b>	Mouse
<b>Immunogen</b>	Human SLAMF7 (aa 249-335).
<b>Isotype</b>	IgG1κ
<b>Purity Detail</b>	Protein G affinity chromatography purified.
<b>Recommendation Dilutions/Conditions</b>	Immunohistochemistry (FFPE) (1-2µg/ml for 30 min at RT)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
<b>Species Reactivity</b>	Human
<b>UniProt ID</b>	Q9NQ25
<b>Worry-free Guarantee</b>	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>

Last modified: July 12, 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)