

S100A12 recombinant monoclonal antibody (S100A12/8952R)

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

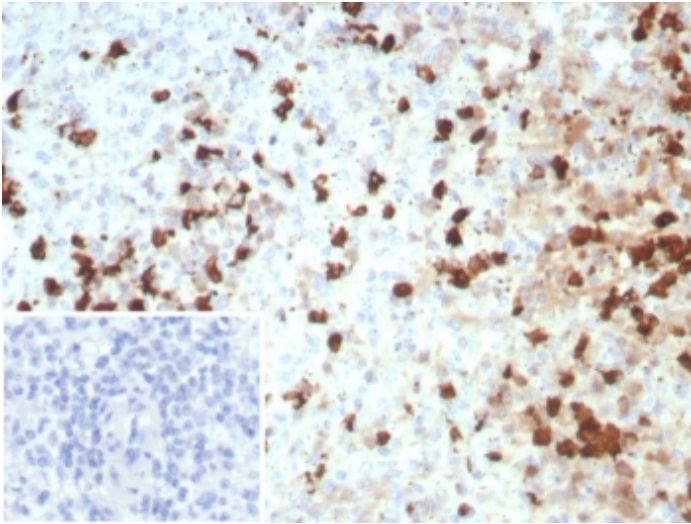
Ordering Information

[Order Online »](#)

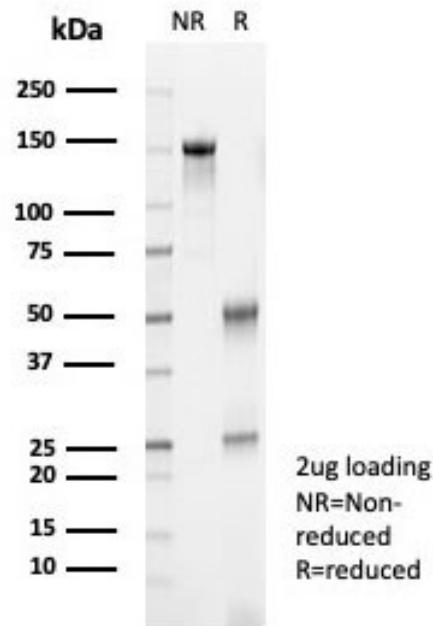
ENZ-ABS918-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



IHC staining of FFPE human spleen tissue with CGRP antibody (clone S100A12/8952R). Inset: PBS used in place of primary Ab (secondary Ab negative control).
 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 S100A12 antibody (clone S100A12/8952R) as confirmation of integrity and purity.

Handling & Storage

Handling	Aliquot the CGRP antibody and store frozen at -20°C or colder. Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	CGRP
Application	IHC (PS)
Clone	S100A12/8952R
Formulation	Liquid. 0.2 mg/ml in 1X PBS containing 0.1 mg/ml BSA (US sourced), 0.05% sodium azide.
Host	Rabbit
Immunogen	A recombinant partial protein sequence (within amino acids 1-95) from human CGRP.
Isotype	IgGκ
Purity Detail	Protein A/G affinity purified.
Recommendation Dilutions/Conditions	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.
Species Reactivity	Human
UniProt ID	P80511
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)