

# S100B Recombinant monoclonal antibody (S100B/1706R)

S100 belongs to the family of calcium binding proteins. S100A and S100B proteins are two members of the S100 family. S100A is composed of an  $\alpha$  and a  $\beta$  chain whereas S100B is composed of two  $\beta$  chains. S-100 protein has been found in normal melanocytes, Langerhans cells, histiocytes, chondrocytes, lipocytes, skeletal and cardiac muscle, Schwann cells, epithelial and myoepithelial cells of the breast, salivary and sweat glands, as well as in glial cells. Neoplasms derived from these cells also express S-100 protein, albeit non-uniformly. A large number of well-differentiated tumors of the salivary gland, adipose and cartilaginous tissue, and Schwann cell-derived tumors express S-100 protein. Almost all malignant melanomas and cases of histiocytosis X are positive for S-100 protein.

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

Citations: 2

[View Online »](#)

## Ordering Information

[Order Online »](#)

|                 |             |
|-----------------|-------------|
| ENZ-ABS307-0100 | 100 $\mu$ g |
| ENZ-ABS307-100  | 100 $\mu$ g |

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

|  |   |
|--|---|
| <b>Alternative Name</b>                    | S100 $\beta$  |
| <b>Application</b>                         | Flow Cytometry, IF, IHC (PS), WB  |
| <b>Application Notes</b>                   | This antibody can be used to localize S-100A and S-100B in various tissue sections.   |
| <b>Clone</b>                               | S100B/1706R   |
| <b>Formulation</b>                         | Liquid. In 1X PBS containing 0.1 mg/ml BSA (US sourced) and 0.05% sodium azide.   |
| <b>Host</b>                                | Rabbit  |
| <b>Immunogen</b>                           | Recombinant full length human S100B.  |
| <b>Isotype</b>                             | IgG $\kappa$  |
| <b>Purity Detail</b>                       | Protein A affinity purified.  |
| <b>Recommendation Dilutions/Conditions</b> | Flow Cytometry (0.5-1 $\mu$ g/10 <sup>6</sup> cells in 0.1 mL)Immunofluorescence (1-2 $\mu$ g/mL)Western blot (0.5-1 $\mu$ g/mL)Immunohistochemistry (formalin-fixed paraffin sections) (0.25-0.5 $\mu$ g/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, Mouse, Rat   |
| <b>Specificity</b>                         | This antibody is specific against an epitope located on the $\beta$ -chain (i.e. in S-100A and S-100B) but not on the alpha-chain of S-100 (i.e. in S-100A and S100A0).   |

UniProt ID

P04271

Worry-free Guarantee

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

Last modified: May 29, 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)