

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 α and a β chain whereas S100B is composed of two β 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 μ g

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name S100 β

Application Flow Cytometry, IF, IHC (PS), WB

Application Notes This antibody can be used to localize S-100A and S-100B in various tissue sections.

Clone S100B/1706R

Formulation Liquid. In 1X PBS containing 0.1 mg/ml BSA (US sourced) and 0.05% sodium azide.

Host Rabbit

Immunogen Recombinant full length human S100B.

Isotype IgGk

Purity Detail Protein A affinity purified.

Recommendation Flow Cytometry (0.5-1 μ g/10⁶ cells in 0.1 mL)Immunofluorescence (1-2 μ g/mL)Western blot (0.5-1 μ g/mL)Immunohistochemistry (formalin-fixed paraffin sections) (0.25-0.5 μ 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.

Dilutions/Conditions

Species Reactivity Human, Mouse, Rat

Specificity This antibody is specific against an epitope located on the β -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

ENZO LIFE SCIENCES

INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES

(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

This antibody is covered by our Worry-Free Guarantee

Belgium, The Netherlands
& Luxembourg

Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France

Phone: +33 1 72 440 1335
info-fr@enzolifesciences.com

Germany

Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland

Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com