

Sp100 (human) polyclonal antibody

Sp100 is an interferon-inducible protein and a major constituent of distinct nuclear domains, known as nuclear dots (NDs) or PML bodies. NDs play a role in autoimmunity, virus infections, and in the etiology of acute promyelocytic leukaemia. SUMO modification of Sp100 has been shown to promote interaction of Sp100 with non-histone chromosomal proteins, suggesting a regulatory role in chromatin organization and in functional interplay between these nuclear bodies and chromatin.

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

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ICC, WB

Formulation Liquid. In PBS containing 0.01% sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to aa 292-304 of human Sp100.

Source From rabbit serum.

Species Reactivity Human

Specificity Recognizes human Sp100 and SUMO modified Sp100 species.

UniProt ID P23497

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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
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