

# SUMO-2 (human) polyclonal antibody

SUMO (small ubiquitin-related modifier) is a member of a family of ubiquitin-like proteins involved in the regulation of the cellular functions of a wide variety of proteins. Four members of the SUMO family have been described in vertebrates, SUMO-1 and the close homologues SUMO-2 and SUMO-3, with some 50% homology between SUMO-1 and SUMO-2/3. Although a fourth SUMO (SUMO-4) has been reported, there remains considerable debate as to whether this is a real and expressed gene product.

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

Citations: 2

[View Online »](#)

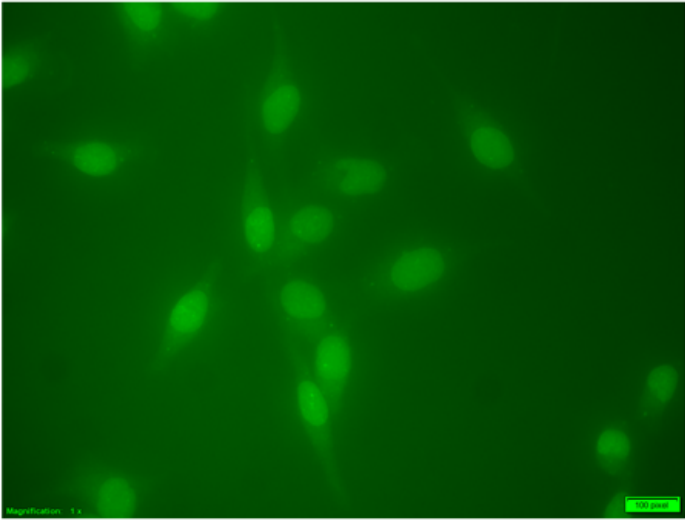
Ordering Information

[Order Online »](#)

|                  |      |
|------------------|------|
| BML-PW0510A-0025 | 25µg |
|------------------|------|

Manuals, SDS & CofA

[View Online »](#)



**Figure 1:** HeLa cells were fixed with paraformaldehyde and stained with SUMO-2 antibody (BML-PW0510A) at 1:50 dilution.

## Handling & Storage

|                           |                                                                                                                                      |
|---------------------------|--------------------------------------------------------------------------------------------------------------------------------------|
| <b>Use/Stability</b>      | Stable for at least 1 year after receipt when stored at -20°C. Store diluted antibody at +4°C, do not freeze and use within 1 month. |
| <b>Handling</b>           | Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.                                                        |
| <b>Short Term Storage</b> | +4°C                                                                                                                                 |
| <b>Long Term Storage</b>  | -20°C                                                                                                                                |
| <b>Shipping</b>           | Blue Ice                                                                                                                             |

**Regulatory Status** RUO - Research Use Only

## Product Details

|                             |                                                                                                                                                  |
|-----------------------------|--------------------------------------------------------------------------------------------------------------------------------------------------|
| <b>Alternative Name</b>     | Sentrin-2, Small ubiquitin-related modifier 2                                                                                                    |
| <b>Application</b>          | ICC, IP, WB                                                                                                                                      |
| <b>Crossreactivity</b>      | Cross-reacts with SUMO-3 due to high sequence homology (86%). Species cross-reactivity expected to be high due to conservation of SUMO sequence. |
| <b>Formulation</b>          | Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.                                                                                   |
| <b>Host</b>                 | Rabbit                                                                                                                                           |
| <b>Immunogen</b>            | Full length recombinant human SUMO-2.                                                                                                            |
| <b>Purity Detail</b>        | Protein A affinity purified.                                                                                                                     |
| <b>Source</b>               | Purified from rabbit serum.                                                                                                                      |
| <b>Species Reactivity</b>   | Human                                                                                                                                            |
| <b>UniProt ID</b>           | P61956                                                                                                                                           |
| <b>Worry-free Guarantee</b> | This antibody is covered by our <a href="#">Worry-Free Guarantee</a>                                                                             |

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)