

BSA (nitrated)

Nitrated bovine serum albumin for use as a positive control for antibodies to nitrotyrosine.

Citations: 1

[View Online »](#)

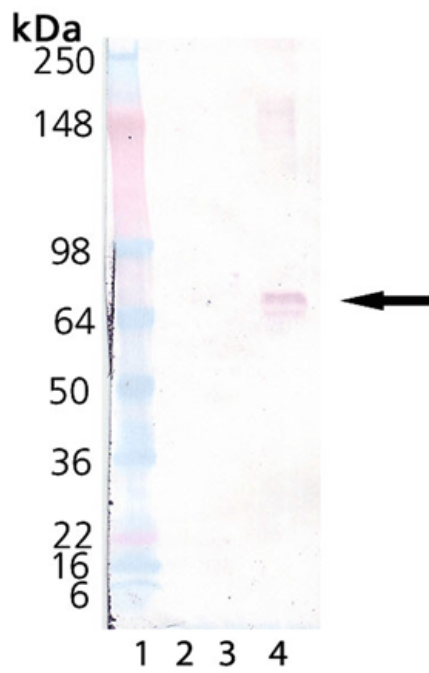
Ordering Information

[Order Online »](#)

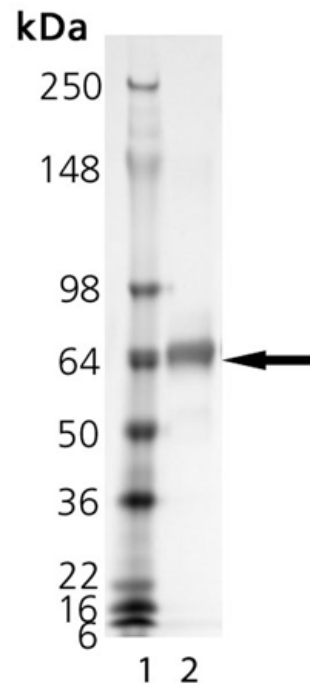
BML-SW125-0100	100µg
----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western Blot analysis: Lane 1: MW Marker, Lane 2: 100 ng BML-SW125, Lane 3: 1 µg untreated BSA, Lane 4: 1 µg BML-SW125. Probed with Nitrotyrosine mAb (Prod no. BML-SE468).



SDS-PAGE Analysis: Lane 1: MW Marker, Lane 2: 1 µg BML-SW125

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Bovine serum albumin

Application Notes For use as a positive control in immunoblots using anti-nitrotyrosine antibodies (Prod. no. BML-SA297, BML-SA468). Use 1-2 µg/lane.

Formulation Lyophilized.

MW ~66kDa

Reconstitution Reconstitute with 200µl of dH2O to yield a 0.5mg/ml solution in PBS.

Source Native. Bovine albumin.

UniProt ID P02769

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-
fr@enzolifesciences.com](mailto:info-fr@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)