

BSA solution (10%), (50 ml)

This BSA Solution (10%) is specially designed to create the Blocking Buffer and Assay Buffer in IMMUNOSET[®] products. Each bottle is sufficient for use with 5x96 wells microtiter plates if following the instructions in our IMMUNOSET[®] protocol. The BSA Solution may be used for any other assay development needs. Additional assay development reagents are available as companion products to our IMMUNOSET[®] products.

Ordering Information

[Order Online »](#)

| | |
|-------------|------|
| ADI-80-1928 | 50ml |
|-------------|------|

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

| | |
|-------------------|--|
| Appearance | Clear to slightly hazy yellowish solution. |
|-------------------|--|

| | |
|--------------------|------------------------------------|
| Formulation | 10% BSA in PBS with preservatives. |
|--------------------|------------------------------------|

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@enzolifesciences.com)

For local distributors and detailed product information visit us online: enzo.com All product names, logos, and brands are the property of their respective owners.