

# BSA solution (10%), (50 ml)

This BSA Solution (10%) is specially designed to create the Blocking Buffer and Assay Buffer in IMMUNOSET<sup>®</sup> products. Each bottle is sufficient for use with 5x96 wells microtiter plates if following the instructions in our IMMUNOSET<sup>®</sup> 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<sup>®</sup> 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.



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)