

# Fixation/Permeabilization (buffer)

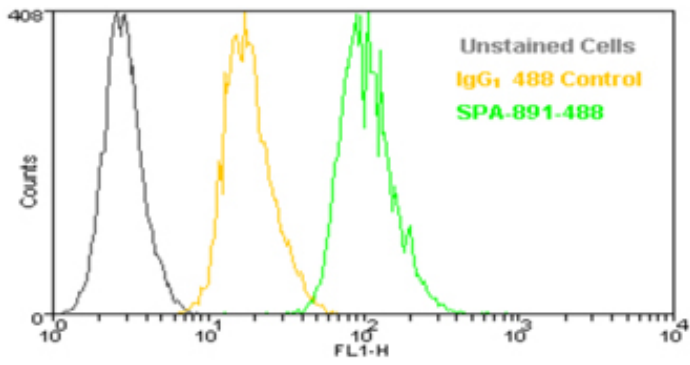
## Ordering Information

[Order Online »](#)

ADI-950-010-100	100ml
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Flow cytometry analysis of 10<sup>6</sup> Jurkat cells were stained using PDI mAb (1D3), DyLight™ 488 Conjugate using one step Fixation/Permeabilization Buffer.

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

### Application Notes

Intended for use in simultaneously fixing and permeabilizing cells in one step to facilitate immunostaining and subsequent analysis of intracellular proteins using flow cytometry. This buffer can be used as an alternative to a two step process with fixation buffer (Prod. No. ADI-950-011) followed by permeabilization buffer (Prod. No. ADI-950-012).

### Formulation

Liquid. PBS, pH 7.2, containing 4% formaldehyde and 0.1% saponin.

### Purity Detail

Purified

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)