

Permeabilization, (buffer)

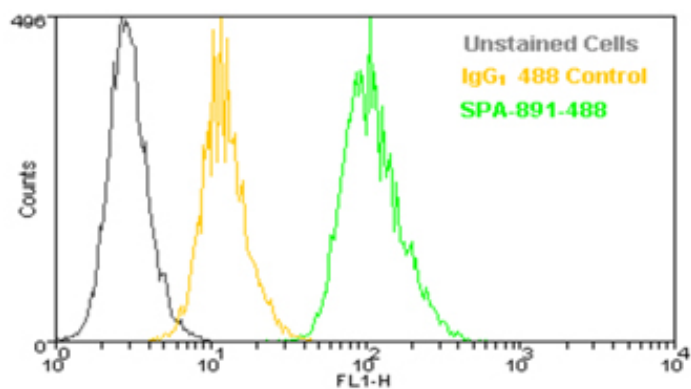
Ordering Information

[Order Online »](#)

ADI-950-012-100	100ml
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Flow cytometry analysis of 10^6 Jurkat cells were stained using PDI mAb (1D3), DyLight™ 488 Conjugate using Fixation Buffer and 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 permeabilizing cells to facilitate immunostaining and subsequent analysis of intracellular proteins using flow cytometry. Use of this buffer should be preceded by fixation of the cells using a suitable fixation buffer (Prod. No. ADI-950-010 or ADI-950-011).

Formulation

Liquid. PBS, pH 7.2, containing 0.1% saponin and 0.09% sodium azide.

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)