

PVDF Transfer Membrane (7cm x 9cm)

Membrane for Western Transfers

PVDF membranes provide high surface area for strong hydrophobic interactions. PVDF membranes typically absorb 50% more protein than nylon or nitrocellulose membranes. This membrane has been optimized for Western transfer applications compatible with WESTERNVIEW® Detection Kits or other chemiluminescent or chromogenic Western blot detection substrates.

Ordering Information

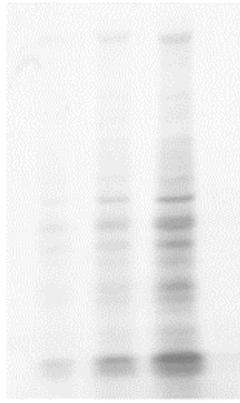
[Order Online »](#)

ENZ-KIT164-0001	10ea
-----------------	------

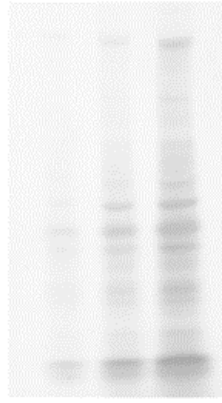
Manuals, SDS & CofA

[View Online »](#)

- Strong protein binding properties
- Sensitive protein detection with low background
- Very low burn-through
- Good chemical compatibility



Enzo
PVDF
Membrane



Competitor
PVDF
Membrane

Rabbit reticulocyte lysate at full strength, 1:3, and 1:10 dilutions. After SDS-PAGE, proteins were transferred to PVDF membranes. Membranes were stained with Amido Black stain for visualization of total protein.

Handling & Storage

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

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)