

PhosphoBlocker™

Blocking Reagent

Most commercially available Western blot blockers, such as dry milk or serum, are sufficient to block unreactive sites on the membrane. However, they are not designed to preserve phosphoprotein antigens during blotting. Our PhosphoBLOCKER™ Blocking Reagent provides superior blocking by maximizing signal-to-noise ratio. The PhosphoBLOCKER™ reagent particularly excels with very low levels of endogenous phosphoproteins.

Ordering Information

[Order Online »](#)

CBI-AKR-104	200g
-------------	------

Manuals, SDS & CofA

[View Online »](#)

- Superior blocking compared to dry milk or serum
- Enhances low level phosphoprotein signal without increasing background
- Preserves phosphoprotein antigens during Western blot procedures
- Easy-to-use premixed dry blend

Handling & Storage

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status

ASR - Analyte Specific Reagent, CE - CE Marked, EUA - Emergency Use Authorization Only, IVD - For In vitro Diagnostic Use Only, IVD - For In vitro Diagnostic Use Only (USA), RUO - Research Use Only, RUO -Research Use Only, Rx Only - For Prescription Use Only

Product Details

Quantity 200 g dry blend; 5% concentration after reconstitution in 4L

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com