



Reveal Your Best Results

Get clear results every time with De-Expose™ background remover for X-ray film.

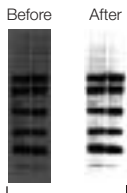
Benefits

- Works on any film, recently exposed or years old
- Doesn't alter results — just gives a cleaner, clearer image
- Requires no reprocessing of gels or blots — treats only the exposed film

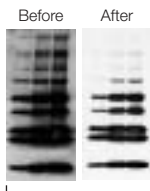


BIO-RAD

Now background can be a thing of the past.



Overexposed film



Overloaded bands

Before and after images demonstrate the effectiveness of De-Expose background remover.

The new De-Expose background remover is available for removal of background from overexposed X-ray film. The kit also removes speckles and corrects overloaded bands. It can be used for any application that uses X-ray film exposure, including western, Southern, and northern blots, gel-shift assays, ribonuclease protection assays, dot blots, and DNA sequencing.

Advantages of De-Expose Background Remover

Never alters results — tested chemistry ensures background is removed from the film, not your data

No prep work required — De-Expose background remover works directly on exposed X-ray film; there is no need to strip, reprobe, and reexpose

Tested for performance on major brands of X-ray film, including Kodak MR, Kodak XAR, and Pierce brands

Works on film even years old so you can retrieve useful data from past projects for current needs

Ordering Information

| Catalog # | Description |
|-----------|--|
| 170-5020 | De-Expose Background Remover, 10x, 250 ml |
| 170-5040 | Immun-Star™ HRP Substrate, 500 ml |
| 170-5041 | Immun-Star HRP Substrate, 100 ml |
| 170-5018 | Immun-Star AP Substrate, 125 ml |
| 162-0115 | Nitrocellulose Membrane, 0.45 μ m, 30 cm x 3.5 m, 1 roll |
| 162-0112 | Nitrocellulose Membrane, 0.2 μ m, 30 cm x 3.5 m, 1 roll |
| 162-0177 | Immun-Blot® PVDF Membrane, 26 cm x 3.3 m, 1 roll |
| 162-0184 | Sequi-Blot™ PVDF Membrane, 26 cm x 3.3 m, 1 roll |
| 162-6404 | Blotting Grade Blocker, nonfat dry milk, 300 g |
| 170-6531 | Tween 20, EIA grade, 100 ml |
| 161-0783 | 1x Phosphate Buffered Saline/1% Casein, 1 L |
| 161-0782 | 1x Tris Buffered Saline/1% Casein, 1 L |

Kodak is a trademark of Eastman Kodak Co. Tween is a trademark of ICI Americas, Inc.

BIO-RAD