



Reveal Your Best Results

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

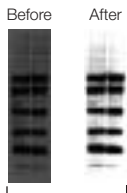
Benefits

- 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

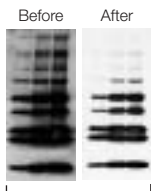


BIO-RAD

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