



# Western blotting at lightning performance.

Western blotting is the most widely used technique for detecting specific proteins. Revvity offers a wide range of western blotting solutions, including our Western Lightning™ and Western Lightning One chemiluminescent substrates. Now, you can have the highest level of sensitivity, specificity, and confidence in your results.

By offering a unique combination of specific immunodetection and size-based separation, western blotting gives you reliable, convenient, high-quality data. Maybe that's why it's still considered the gold standard in protein detection.



## Western Lightning selection guide

Product	Description	Unit size	Part number
<b>Western Lightning ONE - Ready-to-use, premixed single component</b>			
<b>Western Lightning ONE Pico</b>	▪ Mid-picogram detection	25 mL	NEL130001EA
	▪ Ideal for most routine analysis of abundant proteins	250 mL	NEL131001EA
	▪ Low background for high signal to noise ratio	500 mL	NEL132001EA
<b>Western Lightning ONE Femto</b>	▪ Mid-picogram detection	25 mL	NEL130001EA
	▪ Well-suited for a wide range of antibody dilutions	250 mL	NEL131001EA
	▪ Detects protein levels with high signal intensity	500 mL	NEL132001EA
<b>Western Lightning ONE Femto Ultra</b>	▪ Low femtogram detection	25 mL	NEL150001EA
	▪ Low antibody consumption	125 mL	NEL151001EA
	▪ Detects protein levels with superior signal intensity	250 mL	NEL152001EA

Product	Description	Unit size	Part number
<b>Western Lightning - Two-component substrates that require mixing and offer protein detection at varying sensitivities to fit your needs</b>			
<b>Western Lightning Plus</b>	<ul style="list-style-type: none"> <li>▪ Detects proteins in 1-10 picogram range</li> <li>▪ Reduces cost of assay by using fewer primary and secondary antibodies</li> <li>▪ Works well with PVDF and nitrocellulose membranes</li> </ul>	30 mL	NEL103E001EA
		130 mL	NEL103001EA
		340 mL	NEL104001EA
		680 mL	NEL105001EA
<b>Western Lightning Pro</b>	<ul style="list-style-type: none"> <li>▪ Detects proteins in low picogram ranges</li> <li>▪ Delivers outstanding value</li> <li>▪ Offers easy conversion</li> <li>▪ Works well with PVDF and nitrocellulose membranes</li> </ul>	30 mL	NEL120E001EA
		130 mL	NEL120001EA
		340 mL	NEL121001EA
		680 mL	NEL122001EA
<b>Western Lightning Ultra</b>	<ul style="list-style-type: none"> <li>▪ Detects proteins in low femtogram range</li> <li>▪ Uses less of your primary antibody and sample</li> <li>▪ Wide dynamic range, for robust results in just one experiment</li> <li>▪ Immediate, intense signal for at least eight hours; ideal for CCD imagers as well as film</li> <li>▪ Works well with PVDF and nitrocellulose membranes</li> </ul>	20 mL	NEL111001EA
		110 mL	NEL112001EA
		220 mL	NEL113001EA

## Transfer Membranes

Explore Protran™, our high-quality pure nitrocellulose transfer membranes for protein binding. A trusted name in detection, it's easy to use and doesn't require a methanol prewetting step. Our transfer membranes provide superior sensitivity, high resolution, and extremely low background.

Western Blotting Transfer Membranes	Size	Product Description	Part Number
Protran™ Nitrocellulose Transfer Membranes	0.2 µm pore size	30 cm x 3 m Roll	NBA083C001EA
		1 Sheet, 15 cm x 15 cm	NBA083S001EA
		5 Sheets, 33 cm x 56 cm	NBA083G001EA
	0.45 µm pore size	15 cm x 3 m Roll	NBA085A001EA
		20 cm x 3 m Roll	NBA085B001EA
		30 cm x 3.5 m/ Roll	NBA085C001EA
		1 Sheet, 15 cm x 15 cm	NBA085S001EA
		5 Sheets, 33 cm x 56 cm	NBA085G001EA

For more information visit [www.revvy.com](http://www.revvy.com)

