Protein biology

Guide to enzyme substrates for western blotting

Enzyme substrates for western blotting are available in a range of sensitivities for different detection methods. Select a technology and sensitivity to match your experimental requirements.

Chemiluminescent western blotting substrates

Horseradish peroxidase (HRP) substrates

| Thermo Scientific™ substrate | Cat. No. | Signal duration | Detection limit | Primary and secondary antibody dilutions (from 1 mg/mL stock) | Features |
|--|--|--------------------|--------------------------------------|---|--|
| SuperSignal [™] West Atto Ultimate Sensitivity Substrate | A38554 (20 mL) A38555 (100 mL) A38556 (200 mL) | 6 hours | Low femtogram to high attogram | 1°: 1:5,000 2°: 1:100,000 to 1:250,000 | Highest sensitivity substrate for very low-abundance proteins, limited samples, or limited antibodies |
| SuperSignal™ West Femto Maximum Sensitivity Substrate | 34094 (20 mL) 34095 (100 mL) 34096 (200 mL) | 8 hours | Low to mid femtogram | 1°: 1:5,000 2°: 1:100,000 to 1:500,000 | High sensitivity substrate for low- abundance proteins |
| SuperSignal™ West Dura Extended Duration Substrate | 37071 (20 mL) 34075 (100 mL) 34076 (200 mL) | 24 hours | Mid femtogram | 1°: 1:5,000 2°: 1:50,000 to 1:250,000 | Long duration substrate ideal for quantitation |
| SuperSignal [™] West Pico PLUS Chemiluminescent Substrate | 34579 (20 mL) 34577 (200 mL) 34580 (500 mL) 34578 (1 L) | Up to 24 hours | Low picogram to high femtogram | 1°: 1:5,000 2°: 1:20,000 to 1:100,000 | Wide dynamic range for everyday, routine western blotting applications |
| Pierce [™] ECL Plus Western Blotting Substrate | 32134 (25 mL) 32132 (100 mL) | 5 hours | Low picogram to mid femtogram | 1°: 1:1,000 2°: 1:25,000 to 1:200,000 | Can be detected via chemiluminescent and fluorescent imaging technologies; absorption and emission maxima of the substrate are 430 nm and 503 nm, respectively |
| Pierce [™] ECL Western Blotting Substrate | 32109 (50 mL) 32209 (250 mL) 32106 (500 mL) | 1–2 hours | Low picogram | 1°: 1:1,000 2°: 1:5,000 to 1:10,000 | Low cost substrate for high- abundance protein targets |

Alkaline phosphatase (AP) substrates

| Invitrogen [™] substrate | Cat. No. | Recommended membrane | Detection limit | Primary and secondary antibody dilutions (from 1 mg/mL stock) | Maximum light levels reached |
|--|---|-------------------------|----------------------|---|---------------------------------|
| Novex [™] AP Chemiluminescent Substrate (CDP- <i>Star</i> [™]) | WP20002 (100 mL) | PVDF, nylon | Low to mid femtogram | 1°: 1:500 to 1:5,000 2°: 1:2,000 to 1:10,000 | 1 second to 15 minutes |
| CDP- <i>Star</i> [™] Substrate with Nitro-Block-II [™] Enhancer | <u>T2218</u> (100 mL) | Nitrocellulose | Low to mid femtogram | 1°: 1:500 to 1:5,000 2°: 1:2,000 to 1:10,000 | 1 second to 15 minutes |
| CSPD™ Substrate | T2141 (50 mL) T2142 (100 mL) T2143 (250 mL) | PVDF, nylon | High picogram | 1°: 1:500 to 1:5,000 2°: 1:2,000 to 1:10,000 | 2 minutes to 2 hours |
| CSPD™ Substrate with Nitro-Block™ Enhancer | T2217 (100 mL) | Nitrocellulose, PVDF | High picogram | 1°: 1:500 to 5,000 2°: 1:2,000 to 1:10,000 | 2 minutes to 2 hours |



Colorimetric western blotting substrates (precipitating substrates)

HRP substrates

| Thermo Scientific [™] substrate | Cat. No. | Color | Detection limit | Primary and secondary antibody dilutions (from 1 mg/mL stock) | Features |
|--|----------------|-------------|--------------------|---|-------------------------|
| Metal Enhanced DAB Substrate Kit | 34065 (275 mL) | Brown-black | 17 pg | 1°: 1:1,000 2°: 1:5,000 to 1:50,000 | 2-component reagent kit |
| 1-Step™ Ultra TMB Blotting Solution | 37574 (250 mL) | Dark blue | 20 pg | 1°: 1:1,000 2°: 1:5,000 to 1:10,000 | Ready-to-use solution |
| Pierce [™] CN/DAB Substrate Kit | 34000 (275 mL) | Black | 500 pg | 1°: 1:500 2°: 1:2,000 to 1:20,000 | 2-component reagent kit |
| Pierce™ DAB Substrate Kit | 34002 (275 mL) | Brown | 1 ng | 1°: 1:500 2°: 1:2,000 to 1:20,000 | 2-component reagent kit |
| 1-Step™ TMB-Blotting Substrate Solution | 34018 (250 mL) | Dark blue | 1 ng | 1°: 1:500 2°: 1:2,000 to 1:20,000 | Ready-to-use solution |
| 1-Step [™] Chloronaphthol Substrate Solution | 34012 (250 mL) | Blue-purple | 5 ng | 1°: 1:500 2°: 1:2,000 to 1:20,000 | Ready-to-use solution |

AP substrates

| Thermo Scientific [™] substrate | Cat. No. | Color | Detection limit | Primary and secondary antibody dilutions (from 1 mg/mL stock) | Features |
|--|----------------|--------------|--------------------|---|-----------------------|
| 1-Step [™] NBT/BCIP Substrate Solution | 34042 (250 mL) | Black-purple | 30 pg | 1°: 1:500 2°: 1:5,000 to 1:50,000 | Ready-to-use solution |