



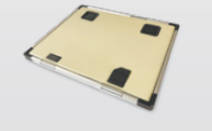
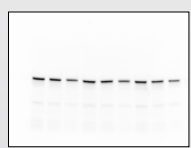
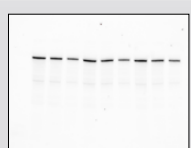
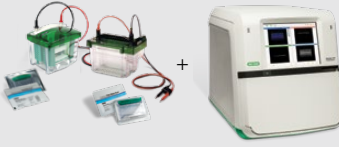
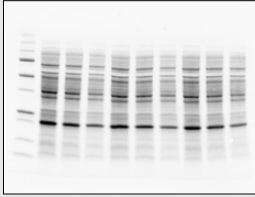

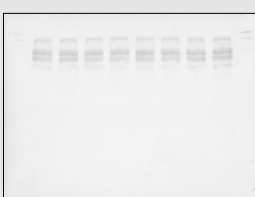

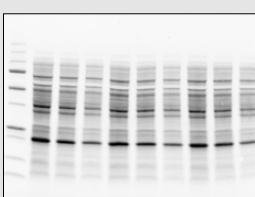
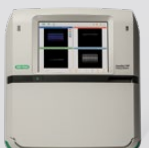
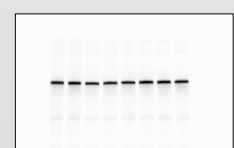
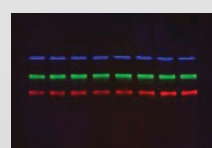


Traditional vs. Stain-Free Western Blotting

Traditional Western Blotting		
Workflow	Time	Data
1 Gel Preparation 	>1 hr gel prep	—
2 Electrophoresis 	~1 hr gel run	—
3 Transfer 	1–3 hr	—
4 Blocking and Antibody Incubation 	~5 hr	—
5 Imaging and Analysis 	>30 min	 Target proteins
6 Strip and Reprobe (Reprobing often needed for actin/tubulin as loading control)	~5 hr	 Loading control
● Total time: 16 hr		

Bio-Rad Stain-Free Western Blotting		
Workflow	Time	Data
1 Electrophoresis TGX Stain-Free Gel, Criterion Cell, ChemiDoc MP Imaging System 	20–30 min	 Pre-transfer gel Stain-Free image to check sample integrity and separation quality
2 Transfer Trans-Blot Turbo Transfer System, ChemiDoc MP Imaging System 	3–10 min	 Post-transfer gel Stain-Free image to measure transfer efficiency
3 Blocking and Antibody Incubation EveryBlot Blocking Buffer, Clarity Western ECL Substrate 	~4 hr	 Stain-Free blot image as loading control
4 Imaging and Analysis ChemiDoc MP Imaging System (No need to strip and reprobe) 	10–15 min	 or  Target proteins (chemiluminescence) Target proteins (fluorescence)
● Total time: 5 hr		

BIO-RAD and TRANS-BLOT are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner.



TGX Stain-Free Precast Gels are covered by U.S. Patent Numbers 7,569,130 and 8,007,646.