




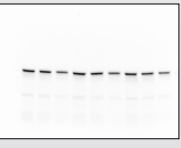
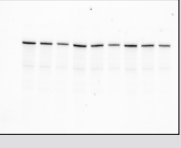
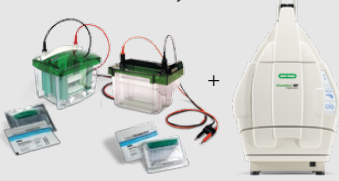
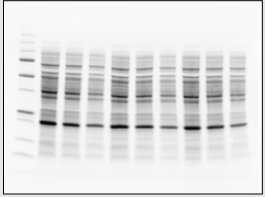

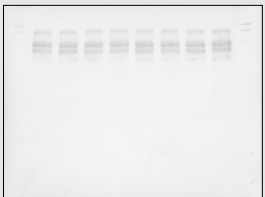

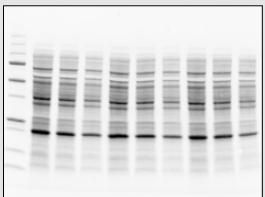

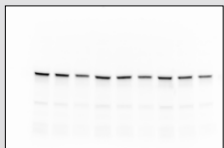
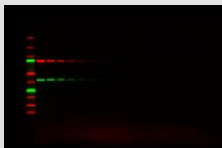


# Traditional vs. V3 Western Workflow™

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

Bio-Rad V3 Western Workflow		
Workflow	Time	Data
<b>1</b> Electrophoresis TGX Stain-Free™ gel, Criterion™ cell, ChemiDoc™ MP system 	20–30 min	 Pre-transfer gel stain-free image to check sample integrity and separation quality
<b>2</b> Transfer Trans-Blot® Turbo™ system, ChemiDoc MP system 	3–10 min	 Post-transfer gel stain-free image to measure transfer efficiency
<b>3</b> Antibody Incubation Clarity™ western ECL substrate 	~5 hr	 Stain-free blot image as loading control
<b>4</b> Imaging and Analysis ChemiDoc MP system (No need to strip and reprobe) 	10–15 min ● Total time: <b>6 hr</b>	 Target proteins (chemiluminescence) or  Target proteins (fluorescence)

