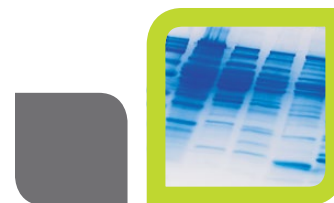
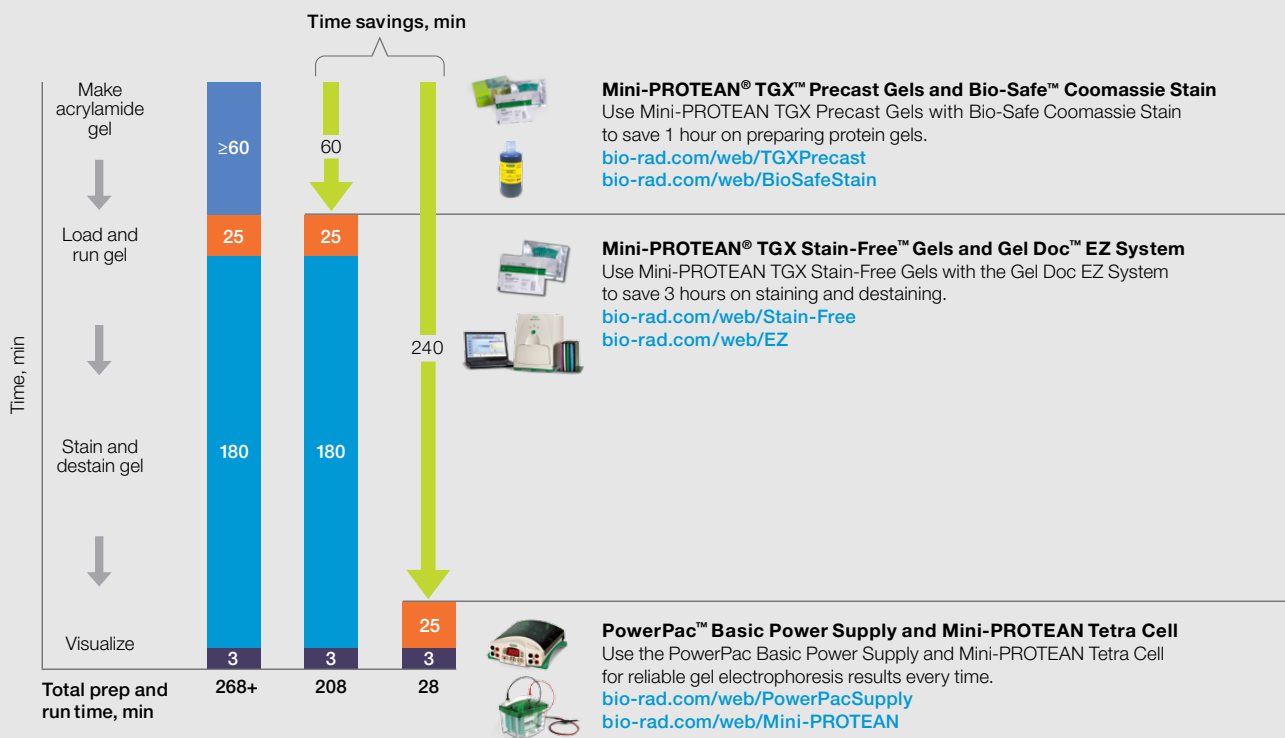


Fast Electrophoresis and Western Blotting

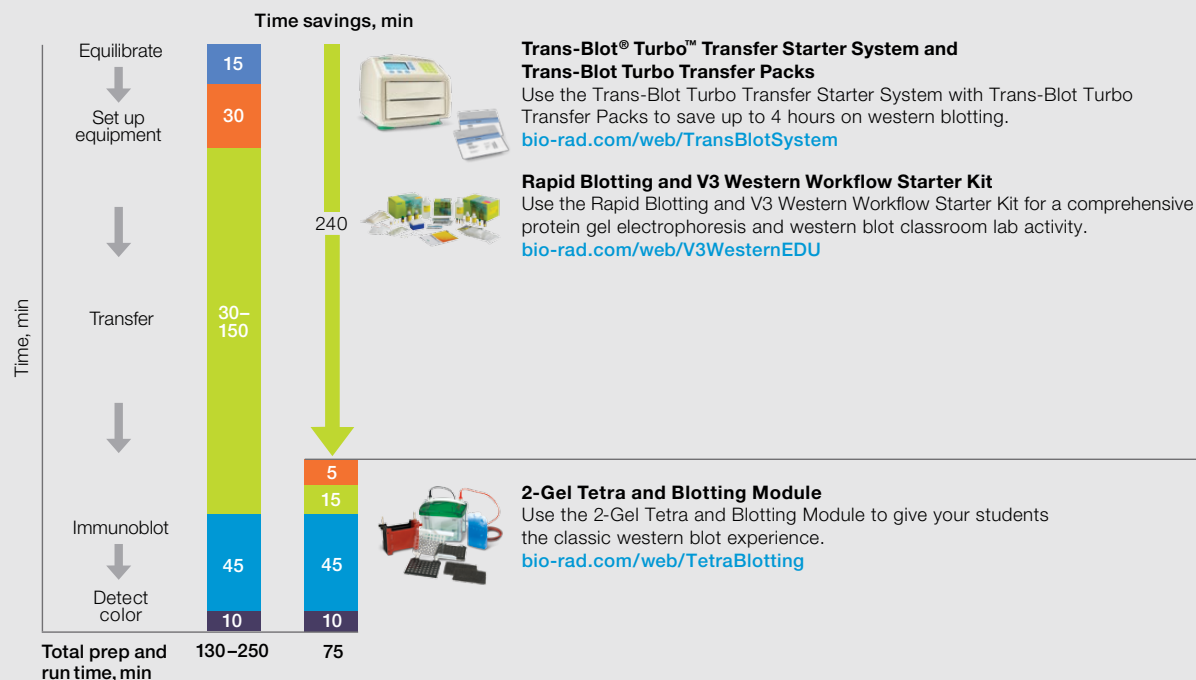


Completing a western blot laboratory in one class can be difficult. But we can help you get more done in less time. Bio-Rad has solutions that save you up to a total of 8 hours — 4 hours with protein gel electrophoresis and 4 hours with western blotting.

Save Up to 4 Hours with Protein Gel Electrophoresis



Save Up to 4 Hours with the V3 Western Workflow™



Visit bio-rad.com/web/V3WorkflowEDU for more information.