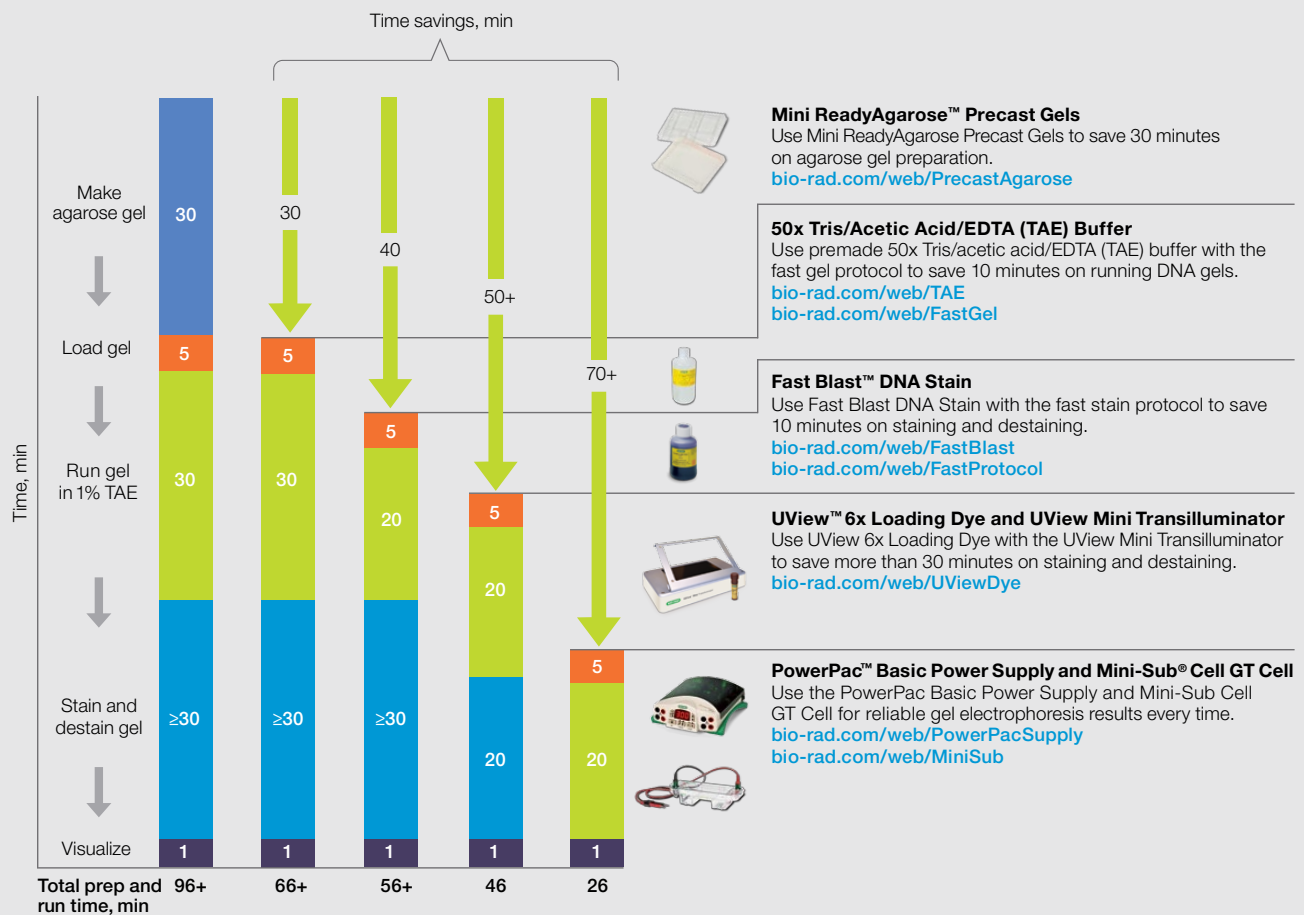


Fast Protein and Nucleic Acid Electrophoresis

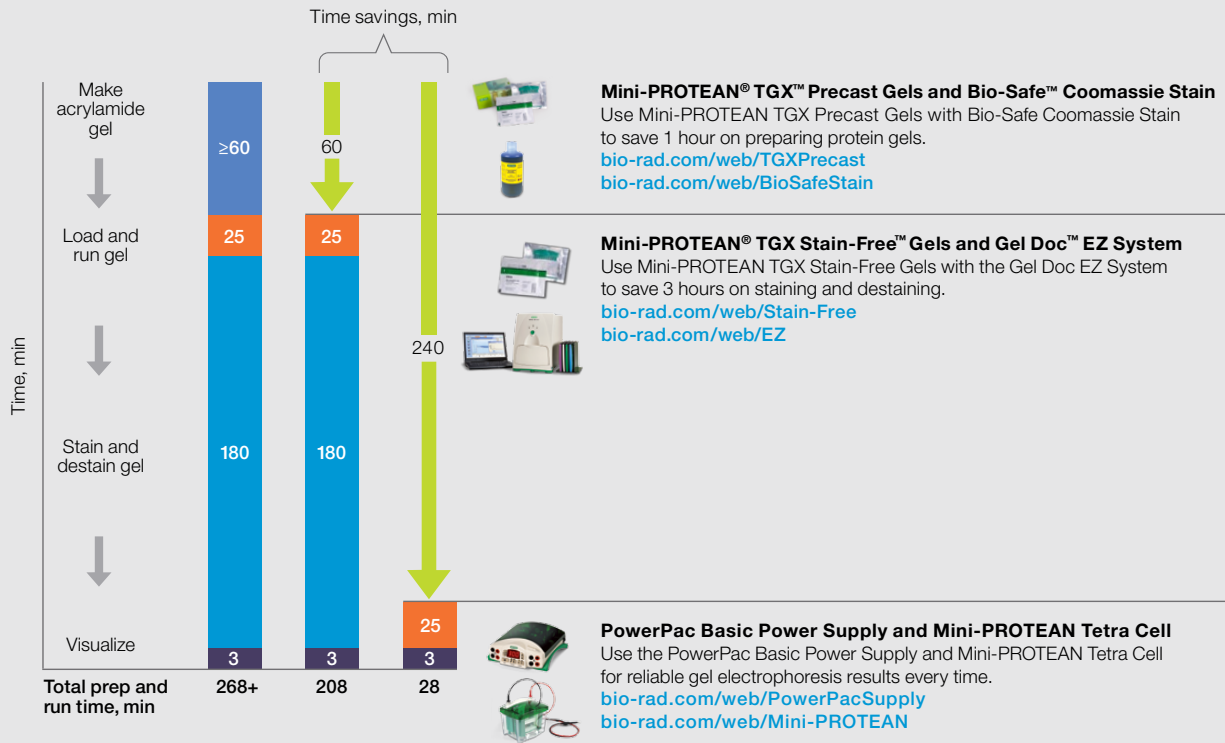


Completing an electrophoresis laboratory in one class can be difficult. We can help you get more done in less time. Bio-Rad has modular solutions to save you more than an hour in preparation and run time for your DNA electrophoresis labs and 4 hours for your protein electrophoresis labs.

Save More than 60 Minutes with DNA Gel Electrophoresis



Save Up to 4 Hours with Protein Gel Electrophoresis



Visit bio-rad.com/web/FastElectroEquipment and bio-rad.com/web/FastElectroReagents for more information about electrophoresis equipment and reagents.