

UView™ 6x Loading Dye and Stain Guide

Use the UView 6x Loading Dye and Stain for all your gel electrophoresis needs!

We've made it easy for you to apply the fast-acting and nontoxic UView Loading Dye and Stain in our line of Biotechnology Explorer™ Kits that use gel electrophoresis. It's easy! Just use the chart below to determine how much UView Loading Dye and Stain to add to the specified kit's DNA marker. There is no need to commit these volumes to memory — UView Loading Dye and Stain is 6x, so just add 1 part to 5 parts sample!

Kit	Catalog Number	DNA Marker Volume, μ l	UView 6x Loading Dye and Stain Added to DNA Marker, μ l	Sample Volume, μ l	UView 6x Loading Dye and Stain Added to Sample, μ l
Cloning and Sequencing Explorer Series	166-5000EDU	400	80	Initial PCR: 20 Nested PCR: 5	Initial PCR: 4 Nested PCR: 1
Fish DNA Barcoding Kit	166-5100EDU	200	40	10	2
PV92 Informatics Kit	166-2100EDU	100	20	Controls: 10 Samples: 20	Controls: 2 Samples: 4
GMO Investigator™ Kit	166-2500EDU	200	40	20	4
Crime Scene Investigator PCR Basics™ Kit	166-2600EDU	200	40	20	4
Forensic DNA Fingerprinting Kit	166-0007EDU	100 (heat marker to 65°C for 5 min, then add UView 6x Loading Dye and Stain)	20	20	4
Restriction Digestion and Analysis of Lambda DNA Kit	166-0002EDU	100 (heat marker to 65°C for 5 min, then add UView 6x Loading Dye and Stain)	20	10	2
Analysis of Precut Lambda DNA Kit	166-0001EDU	100 (heat marker to 65°C for 5 min, then add UView 6x Loading Dye and Stain)	20	10	2