

# Product Information

## 6X DNA Loading Buffer (Blue)

**Catalog Number:** 99962-1

**Unit Size:** 4 x 1.5 mL

### Storage and Handling

Store at 4°C for 6 months or at -20°C for 24 months from date of receipt.

### Product Description

6X Loading Buffer contains density agents for loading DNA samples on agarose or polyacrylamide gels, with two blue electrophoresis tracking dyes that run at approximately 1.5 kb and 200 bp in a 1% agarose gel.

### Protocol

Before gel loading, mix 1 volume of 6X loading buffer with 5 volumes of DNA sample for a final concentration of 1X loading buffer.

Note: The loading buffer does not contain fluorescent DNA dye. For one-step loading and DNA staining, see our 6X GelRed® Prestain Loading Buffers under Related Products. PCR products from EvaGreen® qPCR reactions also can be visualized on gels using a SYBR® Green filter without additional fluorescent dye.

### Related Products

Catalog number	Product Description
31022	Ready-to-Use 1 kb DNA Ladder
31032	Ready-to-Use 100 bp DNA ladder
31039	1 kb DNA Ladder in TE Buffer
31040	100 bp DNA Ladder in TE Buffer
41001	GelRed® Nucleic Acid Gel Stain, 3X in water (4L)
41003	GelRed® Nucleic Acid Gel Stain, 10,000X in water
41005	GelGreen® Nucleic Acid Gel Stain, 10,000X in water
41009	6X GelRed® Prestain Buffer with Blue Tracking Dyes
41010	6X GelRed® Prestain Buffer with Orange Tracking Dye
41008	PAGE GelRed® Nucleic Acid Gel Stain, 10,000X in water
41014	PAGE GelRed® Nucleic Acid Gel Stain, 1X in water (4L)
41020	DNAzure® Visible Blue Nucleic Acid Gel Stain
41006	TBE, 5X (4L)
41024	Water, Ultrapure Molecular Grade (4L)
22020	10X Phosphate Buffered Saline (PBS) (4L)
22007	Activated Charcoal Decontamination Bags
31000	EvaGreen® dye, 20X in water
31041	Forget-Me-Not™ EvaGreen® qPCR Master Mix
31043	Forget-Me-Not™ Universal Probe qPCR Master Mix
31006	AccuBlue® High Sensitivity dsDNA Quantitation Kit
31007	AccuBlue® Broad Range dsDNA Quantitation Kit
31028	AccuClear® Ultra High Sensitivity dsDNA Quantitation Kit
31060	AccuBlue® NextGen dsDNA Quantitation Kit
31066	AccuGreen™ High Sensitivity dsDNA Quantitation Kit for Qubit®
31069	AccuGreen™ Broad Range dsDNA Quantitation Kit for Qubit®

Please visit our website at [www.biotium.com](http://www.biotium.com) for information on our life science research products, including environmentally friendly EvaGreen® qPCR master mixes, DNA quantitation kits, fluorescent CF® dye antibody conjugates and reactive dyes, apoptosis reagents, fluorescent probes, and kits for cell biology research.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.

Qubit and SYBR are registered trademarks of Thermo Fisher Scientific.