

---

# **UView™**

# **6x Loading Dye**

Catalog Numbers

166-5111

166-5112

***BIO-RAD***

## Introduction

Traditionally, a separate dye is used for loading/tracking DNA samples and staining samples for visualization during horizontal electrophoresis. UView™ 6x loading dye simplifies this process because it is a loading dye and fluorescent visualization dye in one formulation. UView dye contains bromphenol blue (comigrates with 500 bp DNA fragments), xylene cyanol (comigrates with 5,000 bp DNA fragments), and a proprietary fluorescent dye for visualization of the DNA after electrophoresis is complete. No staining and destaining are required.

Once the electrophoresis is complete, the gel can immediately be imaged using a UV transilluminator. The fluorescent dye does not interfere with downstream processing of DNA such as restriction digestion or ligation.

Since the fluorescent dye is nontoxic and nonmutagenic, there are no issues with hazardous material shipment fees or hazardous waste disposal fees.

## **Directions for Use**

1. Add 2  $\mu\text{l}$  of UView 6x loading dye to each 10  $\mu\text{l}$  sample of DNA. The final dilution should be 1 part dye to 5 parts DNA sample.
2. Mix the DNA and UView 6x loading dye thoroughly, load samples, and electrophorese. After electrophoresis, place the gel on a UV transilluminator to image the DNA bands. While UView dye enables DNA visualization with ultraviolet light at typical wavelengths of 254, 302, and 365 nm, the optimal wavelength for visualization with UView is 365 nm.

3. UView 6x loading dye is nontoxic and nonmutagenic, so gels can be disposed of according to your state's regulatory requirements.

## Storage Conditions

Store UView 6x loading dye at 4°C.

## Ordering Information

Catalog #	Product Description
161-3101	Certified™ Molecular Biology Agarose, 125 g
166-5111	UView 6x Loading Dye, 200 µl
166-5112	UView 6x Loading Dye, 1 ml
161-0733	TBE Buffer, 10x, 1 Liter
161-0734	TAE Buffer, 50x, 1 Liter
166-0531	UView Mini Transilluminator

***Bio-Rad Laboratories, Inc.***

2000 Alfred Nobel Dr., Hercules, CA 94547 USA