

## Imaging

The 500 bp Fluorescein and Texas Red™ Rulers are easily imaged on the Fluor-S MultiImager system using 2 µl (1 ug) of DNA per lane. When using laser excitation less material (10x less) may be loaded per lane. For multicolor imaging, the 500 bp Fluorescein Ruler can be used with a Texas Red labeled sample and vice versa. DNA labeling kits are available with Fluorescein and Texas Red tagged nucleotides.

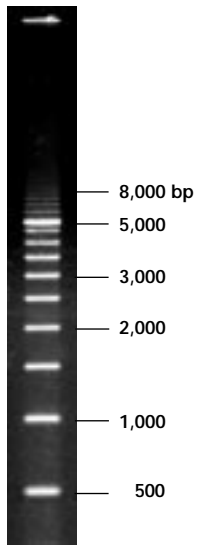


Fig. 1. 1 µg of the 500 bp Texas Red Ruler was diluted to 5 ul in gel loading buffer (2.5% Ficoll, 0.01% bromophenol blue, in TE buffer) and loaded onto a 0.8% agarose gel. The gel was run at 100 V for 90 minutes in 1 X TBE and imaged using 302 nm scanning illumination on the Fluor-S Multilimager system for 2 minutes using the 610 long pass filter.

## 500 bp Texas Red® Ruler

Catalog Number  
**170-8219**

**BIO-RAD**

## Specifications

---

<b>Contents</b>	1 vial 500 bp Texas Red Ruler, 100 ul supplied in TE buffer (10 mM Tris-HCl, 1 mM EDTA, pH 8). 1 vial 5X Gel Loading Buffer (5X TE, 15% Ficoll, 0.05% Bromophenol Blue).
<b>Quantity</b>	DNA sufficient for 50 lanes when used at 2 ul per lane.
<b>Concentration</b>	500 ug/ml
<b>Shipping</b>	The 500 bp Texas Red Ruler is shipped at room temperature.
<b>DNA fragment sizes</b>	The 500 bp Texas Red Ruler contains 16 bands, 500 to 8,000 bp in exact 500 bp increments. A visibly brighter reference band is present at 5,000 bp.

## Storage

The 500 bp Texas Red Ruler should be stored at 4 °C in the dark. For long term storage the standard can be stored at -20 °C Use only sterile pipette tips when removing aliquots. Introduction of nucleases will shorten the shelf life.

## Shelf life

The 500 bp Texas Red Ruler is stable for 1 year when stored in the dark at 4 °C.

## Optimum conditions

The 500 bp ruler can be well resolved on a 0.4% to 2% agarose gel when the appropriate gel length is considered. Most agarose has a fluorescent component and using Bio-Rad

Molecular Biology Certified Agarose will minimize the background effects associated with the fluorescence in agarose. Also, a thin gel (3 mm) will have less background than a thick gel (> 5 mm).

Example, an 0.8% agarose gel, 3mm thick and 10 cm in length, should be run at 90 to 100 volts for 1.5 hours or until the dye marker is at least 3/4 of the way down the gel. This will allow the dye marker to move past the 500 bp band, and allow for good resolution of the bands.