

passes light in the range of 600 to 650 nm on other imaging devices. For multicolor imaging, we recommend labeling of the unknown with fluorescein and electrophoresing an equivalent amount of the Texas Red-labeled standard. Virtually no signal is seen from Texas Red when imaging fluorescein with the 530DF60 band pass filter on the Fluor-S MultiImager system .

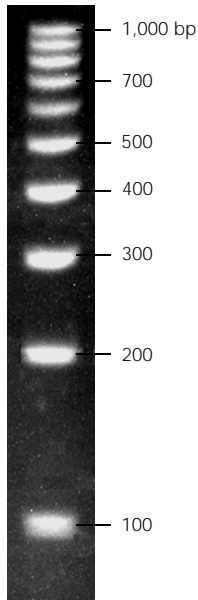


Fig. 1. 400 ng of the 100 bp Texas Red Ruler was diluted to 5 ul in gel loading buffer (2.5% Ficoll, 0.01% bromophenol blue, in TE buffer) and loaded onto a 5% polyacrylamide gel. The gel was run at 100 V for 60 minutes in 1x TBE and imaged using 302 nm scanning illumination on the Fluor-S MultiImager for 30 seconds with the 610 long pass filter.



100 bp Texas Red® Ruler

**Catalog Number
170-8217**



Specifications

Contents	1 vial 100 bp Texas Red Ruler, 100 µl 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).
Quantity	DNA sufficient for 50 lanes when used at 2 µl per lane.
Concentration	200 µg/ml
Shipping	The 100 bp Texas Red Ruler is shipped at room temperature.
DNA fragment sizes	The 100 bp Texas Red Ruler contains 10 visible bands, 100 to 1,000 bp in exact 100 bp increments.

Storage

The 100 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 100 bp Texas Red Ruler is stable for 1 year when stored in the dark at 4 °C.

Use

The 100 bp Texas Red Ruler can be well resolved on polyacrylamide gels of less than 10% or on high concentration, > 2.5%, agarose gels. For imaging with 302 nm (UV) excitation, *e.g.* the Bio-Rad Fluor-S™ MultiImager system, 2 µl (400 ng) of the DNA should be loaded per lane, following dilution in gel loading buffer. When using laser excitation, less material (~10 times less) may be loaded per lane. Adjustments may be made to the loading volume for different well sizes and desired band intensity. The Texas Red label is best imaged with the 610 long pass filter on the Fluor-S MultiImager system or with a filter which optimally