

---

## trUView™ Cuvettes

Disposable UV Transparent Cuvettes

### Catalog Numbers

170-2510

170-2511

170-2512

### Description

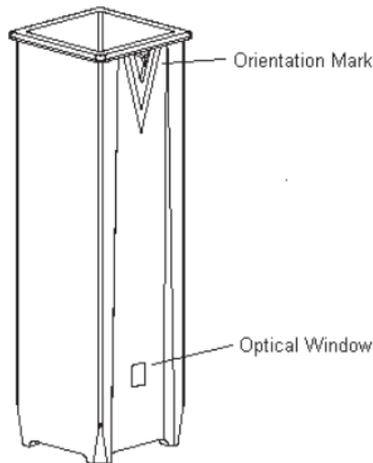
trUView Cuvettes, 50

trUView Cuvettes, 100

Height Adaptor, 15 mm

## Introduction

The trUView™ cuvette can supplement or replace the use of quartz cuvettes for the quantitation of common molecular biology samples. Ample light transmission in the UV wavelengths provides for accurate quantitation of nucleic acid samples.



## Principle of Operation

The trUView cuvette is a disposable plastic cuvette suitable for molecular biology assays from 220 to 1,100 nm. The trUView cuvette is constructed from plastic that is highly transmissible in both the UV and visible light range. While

many other types of plastic absorb light at wavelengths below 400 nm, the trUView cuvette maintains excellent transmission of light at 260 nm, permitting quantitation of nucleic acids.

The trUView cuvette is compatible with most spectrophotometers and no height adaptor is required for those instruments with a light path height (z-dimension) of 8.5 mm. A height adaptor is available from Bio-Rad for use with spectrophotometers with light path heights of 15 mm. Consult your spectrophotometer's user manual for z-dimension information.

Use the trUView cuvette to completely replace quartz cuvettes or as a supplement to their use when working with precious DNA or RNA samples. These cuvettes are certified DNase and RNase free, so the sample may be recovered from the cuvette and used for additional assays. Sample volumes as low as 50 µl may be used.

## Operating Instructions

trUView cuvettes are compatible with spectrophotometers having a light-path height (z-dimension) of 8.5 mm. Consult your spectrophotometer's user manual for z-dimension information.

For most spectrophotometers, use the following protocol to ensure the most accurate assay results.

1. Ensure that the two optical windows of the cuvette are clean and free from scratches.
2. **Determine the baseline absorbance (blank) using a trUView cuvette.** For the most accurate results, be sure to note the orientation of the cuvette (marked with an arrow at the top of the cuvette) and use a buffer that is identical to that used with the sample.
3. Insert the trUView cuvette with 50 µl of sample into the cuvette chamber in the same orientation as used when determining baseline absorbance.

## Other Manufacturers

Instruments with a light path height of 15 mm will require the use of an optional adaptor (170-2512). The height adaptor is inserted into the cell holder of the spectrophotometer prior to the assay. After the adaptor has been inserted, the standard protocol may be followed.

## Troubleshooting

### Inaccurate Assay Values

- Ensure blank and sample are assayed using the same cuvette orientation.
- For best results measure the blank and sample in the same trUView cuvette.
- Inspect the cuvette for scratches or dirt on optical windows.
- Ensure the light path height (z-dimension) of the spectrophotometer is 8.5 mm. If it is 15 mm, use the optional height adaptor.
- If cuvette has been reused, discard and use a new cuvette. Multiple assays will degrade light transmission properties of the cuvette.
- Visually check the sample has no air bubbles. Air bubbles can be removed by gently tapping the cuvette.

## Technical Specifications

Outer Dimensions	12.5 x 12.5 x 45 mm
Light Path Height	8.5 mm
Pathlength	10 mm
Functional Wavelength Range	220–1,100 nm
Baseline Absorbance	$\leq 0.3$ AU @ 260 nm
Variability	$\leq 15$ mAU @ 260 nm
Functional Volume	$\geq 50$ $\mu$ l

### Chemical Compatibility

trUView cuvettes are resistant to all chemical buffers typically used in most life science laboratories. The trUView cuvettes are not compatible with some solvents such as chloroform and toluene.

### Contaminants

All trUView cuvettes are guaranteed RNase and DNase free.