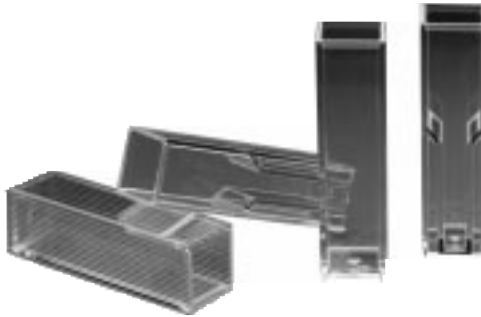


# Disposable Cuvettes

**Visible Range, 800-340 nm  
Standard and Semi-micro,  
for use with Bio-Rad  
Protein Assay Kits**



For Technical Service  
Call Your Local Bio-Rad Office or  
in the U.S. Call **1-800-4BIORAD**  
(1-800-424-6723)

**BIO-RAD**



# Introduction

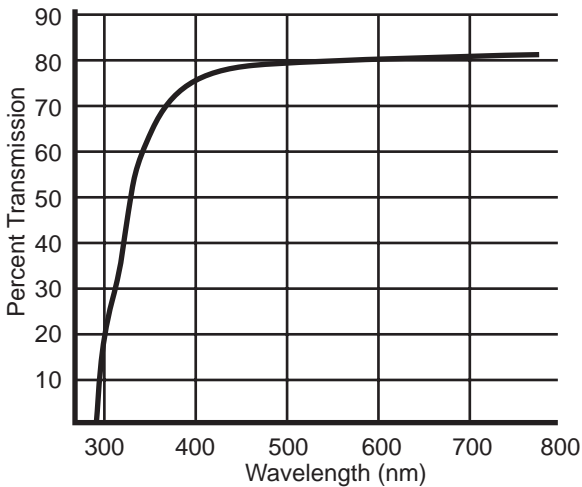
Why use expensive glass or quartz cuvettes for many applications, such as protein assay, where you can use excellent quality, disposable polystyrene cuvettes? Bio-Rad's cuvettes fit virtually all spectrophotometer standard cuvette holders, including Beckman®, Gilford®, Hatachi, Perkin-Elmer, Varian, and Zeiss.

- Excellent specifications
- Fit most spectrophotometers
- Delivery from stock
- Inexpensive
- Unconditionally guaranteed by Bio-Rad

## Specifications

---

<b>Energy levels</b>	≥ 80% T at 800 nm ≥ 55% T at 340 nm
<b>Transmission matching</b>	± 1%
<b>Path length matching</b>	± 0.5%
<b>Path length</b>	1 cm
<b>Material</b>	Premium grade virgin polystyrene
<b>Dimensions</b>	12.5 mm square x 45 mm high
<b>Capacity</b>	
<b>Standard</b>	3.5 ml
<b>Semi-micro</b>	1.5 ml



The performance specifications were obtained using a Perkin-Elmer UV/VIS spectrophotometer (scanning double-beam instrument plus output device) with empty cuvettes versus air path and no modifications to 'improve' results.

## Disposable Cuvettes for Protein Assays

Bio-Rad's disposable cuvettes adsorb less Coomassie blue dye than glass cuvettes, and much less than quartz. We recommend the standard cuvettes for use with the Bio-Rad Protein Assay (based on the Bradford method) and the Bio-Rad *DC* (detergent-compatible) Protein Assay Kits. The cuvettes can be used as mixing vessels, eliminating messy transfer and clean-up steps between sample measurements. In addition, the volume of reagents can be cut in half, yielding twice as many assays per kit. (For additional information on protein assays, request bulletin 1069.)

The low cost of matched cuvettes allows the use of one cuvette for each sample. This eliminates time consuming, repetitive transfer of samples to spectrophotometer cells.

While optical surfaces are precisely smooth to yield the performance indicated, non-optical walls are grooved from top to bottom to provide an easy grip and quick orientation, and to discourage accidental fingerprinting of optical surfaces or misorientation in machines.



Cuvettes are packaged 100 per tray. Each dust proof styrofoam tray has its own lid and is packed in a locking carton for shipment. In this way you never need expose even one tray of cuvettes to ambient dust, as happens when one styrofoam block serves as both tray and lid in a 'sandwich' arrangement.

# Ordering Information

<b>Catalog Number</b>	<b>Product Description</b>
223-9950	<b>Disposable Polystyrene Cuvettes Standard</b> , 3.5 ml, 100/covered, compartmented styrofoam tray
223-9955	<b>Disposable Polystyrene Cuvettes Semi-micro</b> , 1.5 ml, 100/covered, compartmented styrofoam tray

**BIO-RAD**

**Bio-Rad  
Laboratories**

*Life Science  
Group*

**U.S.** (800) 4BIORAD • **California** Ph. (510) 741-1000 • **New York**  
Ph. (516) 756-2575 • **Australia** Ph. 02-805-5000 • **Austria**  
Ph. (1) 877 89 01 • **Belgium** Ph. 09-385 55 11 • **Canada**  
Ph. (905) 624-0713 • **China** Ph. (01) 2563148 • **France** Ph. (1) 49 60 68 34  
• **Germany** Ph. 089 31884-0 • **Italy** Ph. 02-21609 1 • **Japan**  
Ph. 03-3534-7515 • **Hong Kong** Ph. 7893300 • **The Netherlands**  
Ph. 08385-40666 • **New Zealand** Ph. 09-443 3099 • **Singapore**  
Ph. (65) 4432529 • **Spain** Ph. (91) 661 70 85 • **Sweden**  
Ph. 46 (0) 735 83 00 • **Switzerland** Ph. 01-810 16 77 • **United  
Kingdom** Ph. 0800 181134

SIG 093093 Printed in USA

LIT254 Rev C