

Liquid Handling

Pipets and Disposables for Microplate Systems



Products	Features and Benefits
Costar 96-well flat-bottom EIA plates	<ul style="list-style-type: none"> ▪ High binding ▪ High optical clarity ▪ Low background absorption
Sterilized reagent reservoirs	<ul style="list-style-type: none"> ▪ Ideal for dispensing reagent with multichannel pipets ▪ Sloping design makes it easy to pick up liquid to the last drop
Micro test tubes	<ul style="list-style-type: none"> ▪ Suitable for general-purpose benchtop centrifuge use ▪ Autoclavable to 120°C; freezable to -80°C
Xcluda™ aerosol barrier pipet tips	<ul style="list-style-type: none"> ▪ Guard against aerosol contamination of samples ▪ Certified to be free of DNase, RNase, and pyrogens
Standard pipet tips	<ul style="list-style-type: none"> ▪ Accurately molded for an airtight fit ▪ All materials formulated without heavy metals



Ordering Information

Catalog # Description

96-Well EIA Plate

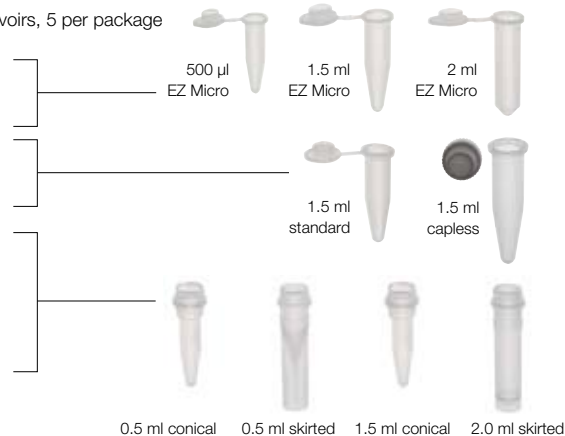
224-0096 **Costar 96-Well Flat-Bottom EIA Plates**, polystyrene, box of 100 plates, 5 per package

Reagent Reservoirs

224-4872 **Sterilized Reagent Reservoirs**, graduated, polystyrene, box of 200 reservoirs, 5 per package

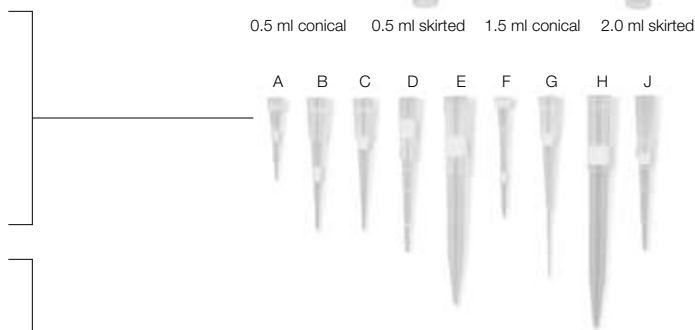
Micro Test Tubes

- 223-9503 **EZ Micro™ Test Tubes**, 500 µl, natural, 1,000
- 223-9480 **EZ Micro Test Tubes**, 1.5 ml, natural, 500
- 223-9430 **EZ Micro Test Tubes**, 2 ml,* natural, 500
- 223-9501 **Micro Test Tubes**, standard, 1.5 ml, natural, 500
- 223-9500 **Micro Test Tubes**, capless, 1.5 ml, polypropylene, natural, graduated, 500
- 223-9490 **Separate Caps**, for capless micro test tubes, white, 1,000
- 224-0165 **0.5 ml Conical Tubes**, with installed O-ring screwcaps, sterilized, 500
- 224-0185 **0.5 ml Skirted Tubes**, with installed O-ring screwcaps, sterilized, 500
- 224-0100 **1.5 ml Conical Tubes**, with separate O-ring screwcaps, unsterilized, 500
- 224-0110 **1.5 ml Conical Tubes**, with installed O-ring screwcaps, sterilized, 500
- 224-0130 **2.0 ml Skirted Tubes**, with separate O-ring screwcaps, unsterilized, 500
- 224-0140 **2.0 ml Skirted Tubes**, with installed O-ring screwcaps, sterilized, 500



Xcluda Aerosol Barrier Pipet Tips

- 211-2001 **Xcluda Style A**, 0.5–10 µl, 960
- 211-2006 **Xcluda Style B**, 2–20 µl, 960
- 211-2011 **Xcluda Style C**, 10–100 µl, 960
- 211-2016 **Xcluda Style D**, 20–200 µl, 960
- 211-2021 **Xcluda Style E**, 100–1,000 µl, 960
- 211-2026 **Xcluda Style F**, 0.1–10 µl, 1,000
- 211-2031 **Xcluda Style G**, 10–100 µl, 1,000
- 211-2036 **Xcluda Style H**, 100–1,000 µl, 1,000
- 211-2041 **Xcluda Style J**, 10–100 µl, 960



Standard Pipet Tips

- 223-9014 **BR-14 Tips**, natural, 0.1–10 µl, 1,000
- 223-9026 **BR-26 Tips**, natural, 5–200 µl, 1,000
- 223-9028 **BR-28 Tips**, natural, 25–200 µl, 1,000
- 223-9031 **BR-31 Tips**, natural, 0.5–10 µl, 1,000
- 223-9035 **BR-35 Tips**, yellow, 1–200 µl, 1,000
- 223-9037 **BR-37 Tips**, natural, 1–200 µl, 1,000
- 223-9038 **BR-38 Tips**, yellow, graduated, beveled, 1–200 µl, 1,000
- 223-9039 **BR-39 Tips**, natural, graduated, beveled, 1–200 µl, 1,000
- 223-9040 **BR-40 Tips**, blue, 100–1,000 µl, 500
- 223-9041 **BR-41 Tips**, natural, 100–1,000 µl, 500



Racked Pipet Tips, One Cover on Nested Racks

223-9135 **RBR-35 Tips**, yellow, 1–200 µl, 1,000

Racked Pipet Tips, Cover on Each Rack

- 223-9354 **TBR-14 Tips**, natural, 0.1–10 µl, 1,000
- 223-9347 **TBR-35 Tips**, yellow, 1–200 µl, 1,000
- 223-9350 **TBR-40 Tips**, blue, 100–1,000 µl, 1,000
- 223-9351 **TBR-41 Tips**, natural, 100–1,000 µl, 1,000

* To fit centrifuges, 2 ml tubes have thinner walls than standard tubes and are intended for short runs at less than 13,000 x g. Heavy samples, high g-forces, long runs, or centrifuges that radiate excess heat may damage these tubes.

For more information, go to www.bio-rad.com/liquidhandling/

Costar is a trademark of Corning Costar Corporation.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA 800 4BIORAD Australia 61 02 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3237 9400
Canada 905 364 3435 China 86 21 6426 0808 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65
Germany 089 318 84 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 455 8800 India 91 124 4029300 Israel 03 963 6050
Italy 39 02 216091 Japan 03 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666 New Zealand 0508 805 500
Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 495 721 14 04 Singapore 65 6415 3188 South Africa 27 861 246 723
Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan 886 2 2578 7189 United Kingdom 020 8328 2000