



MATTEK 
A BICO COMPANY

PRODUCT CATALOG

MatTek Cultureware





Summary

About MatTek

MatTek Dishes

Dish Properties

Glass Bottom Multi-Well Plates

Coverslips and Coverslip Kits

Chambered Cell Culture Slides

PermaCell Cell Culture Inserts

Cultureware FAQ's

Supporting Your Science

Founded in 1985, MatTek Life Sciences was created by two MIT chemical engineering professors as a cell culture surface technology company. In 1993, MatTek began producing its first lab-grown human skin tissue model, EpiDerm, as a reliable replacement for animal testing. MatTek has gone on to produce additional human organ models including EpiOcular, EpiOral, EpiGingival, EpiVaginal, EpiAirway, EpiAlveolar, EpiNasal, and EpiIntestinal. Since 2004, EpiDerm has achieved worldwide regulatory acceptance and validation validations for Skin Irritation (OECD TG 439), Skin Corrosion (OECD TG 431), Phototoxicity (OECD TG 498), and Skin Irritation for Medical Device Extracts (ISO10993:2021). These validations not only signified the reliability and relevance of EpiDerm for preclinical testing but also the broad gap that exists between animal models and human biology. In addition to steady growth in their U.S. location, MatTek established a European location in Bratislava, Slovakia, in 2009 to expand the global availability of its human tissue technology.

In March 2021, MatTek was acquired by BICO (formerly CELLINK), the leading global bio-convergence company with headquarters in Gothenburg, Sweden. This extraordinary group of scientists shares MatTek's vision for the future of our industry and technologies. Today, MatTek's physiologically advanced 3D tissue models of the skin, eye, oral, respiratory, vaginal, and intestinal systems empower companies in the cosmetics, chemical, pharmaceutical, and consumer goods industries to assess the safety and efficacy of their formulations, chemicals, and compounds without the use of animals. MatTek is proud to offer non-animal testing platforms that lower preclinical costs and provide more human-relevant results. In addition to its human tissue technology, MatTek offers a selection of primary cells, customizable culture media, 3DIY Tissue Kits, in-house testing services, and glass bottom MatTek Dishes for brilliant microscopic imaging. We are here to support your science every step of the way.

MatTek Dishes

MatTek Glass Bottom Dishes combine the convenience of standard size 35 mm, 50 mm, 60 mm, and 100 mm plastic Petri dishes with the optical quality of glass, providing superior microscopic images. A removable coverslip allows for increased imaging flexibility.



Optional coatings including poly-D-lysine and type 1 rat tail collagen that assist with cell adherence to the glass are available as standard products.

STERILE

All MatTek dishes are sterilized with gamma-irradiation

GLASS DIAMETER

| | 7 mm | 10 mm | 14 mm | 20 mm | 30 mm |
|--------|------|-------|-------|-------|-------|
| 35 mm | ■ | ■ | ■ | ■ | |
| 50 mm | | | ■ | | ■ |
| 60 mm | | | | ■ | ■ |
| 100 mm | | | | | ■ |

For most applications, a 14 mm glass microwell is sufficient. Choose the 20 mm diameter microwell to maximize viewing area or a 10 mm diameter if your application is cell-, media-, or reagent-limited.

COVERSLIP

| | No. 0 | No. 1.0 | No. 1.5 | No. 0.170 |
|--------|-------|---------|---------|-----------|
| 35 mm | ■ | ■ | ■ | ■ |
| 50 mm | ■ | | ■ | |
| 60 mm | | | ■ | |
| 100 mm | | | ■ | |

For most applications, No. 1.5 is the preferred coverslip thickness, especially for optimizing image quality with high numerical aperture objectives. The No. 0 coverslip gives you the most working distance and may be useful for thicker specimens.



35 mm MatTek Dishes

Everyone's favorite dish for microscopy! MatTek's 35 mm Glass Bottom Dishes combine the convenience of standard size, disposable plastic Petri dishes with the optical quality of glass, providing superior microscopic images.

PART NO.

DESCRIPTION

35 mm Dish | No. 0 Coverslip

10 mm Glass Diameter

| | |
|----------------|--|
| P35G-0-10-C | No. 0 Coverslip, 10 mm Glass Diameter |
| P35GC-0-10-C | No. 0 Poly-D-Lysine Coated Coverslip, 10 mm Glass Diameter |
| P35GCOL-0-10-C | No. 0 Collagen Coated Coverslip, 10 mm Glass Diameter |

14 mm Glass Diameter

| | |
|----------------|--|
| P35G-0-14-C | No. 0 Coverslip, 14 mm Glass Diameter |
| P35GC-0-14-C | No. 0 Poly-D-Lysine Coated Coverslip, 14 mm Glass Diameter |
| P35GCOL-0-14-C | No. 0 Collagen Coated Coverslip, 14 mm Glass Diameter |

7 mm and 20 mm Glass Diameter

| | |
|-------------|--------------------------------------|
| P35G-0-7-C | No. 0 Coverslip, 7 mm Glass Diameter |
| P35G-0-20-C | No. 0 Coverslip, 20mm Glass Diameter |

35 mm Dish | No. 1.0 Coverslip

14 mm Glass Diameter

| | |
|------------------|--|
| P35G-1.0-14-C | No. 1.0 Coverslip, 14 mm Glass Diameter |
| P35GC-1.0-14-C | No. 1.0 Poly-D-Lysine Coated Coverslip, 14 mm Glass Diameter |
| P35GCOL-1.0-14-C | No. 1.0 Collagen Coated Coverslip, 14 mm Glass Diameter |

20 mm Glass Diameter

| | |
|---------------|---|
| P35G-1.0-20-C | No. 1.0 Coverslip, 20 mm Glass Diameter |
|---------------|---|

35 mm Dish | No. 1.5 Coverslip

7 mm Glass Diameter

| | |
|--------------|--|
| P35G-1.5-7-C | No. 1.5 Coverslip, 7 mm Glass Diameter |
|--------------|--|

10 mm Glass Diameter

| | |
|------------------|--|
| P35G-1.5-10-C | No. 1.5 Coverslip, 10 mm Glass Diameter |
| P35GC-1.5-10-C | No. 1.5 Poly-D-Lysine Coated Coverslip, 10 mm Glass Diameter |
| P35GCOL-1.5-10-C | No. 1.5 Collagen Coated Coverslip, 10 mm Glass Diameter |

14 mm Glass Diameter

| | |
|------------------|--|
| P35G-1.5-14-C | No. 1.5 Coverslip, 14 mm Glass Diameter |
| P35GC-1.5-14-C | No. 1.5 Poly-D-Lysine Coated Coverslip, 14 mm Glass Diameter |
| P35GCOL-1.5-14-C | No. 1.5 Collagen Coated Coverslip, 14 mm Glass Diameter |

20 mm Glass Diameter

| | |
|---------------|---|
| P35G-1.5-20-C | No. 1.5 Coverslip, 20 mm Glass Diameter |
|---------------|---|

35 mm Dish | No. 1.5 Coverslip (High Precision & Additional)

14 mm Glass Diameter

| | |
|--------------------|--|
| P35G-0.170-14-C | No. 1.5 (High Precision) Coverslip, 14 mm Glass Diameter |
| P35G-1.5-14-C-HA | No. 1.5 Coverslip, 14 mm Glass Diameter, Treated for High Adhesion |
| P35P-1.5-14-C | No. 1.5 Plastic Coverslip, 14 mm Plastic Diameter |
| P35G-1.5-14-C-GRD | No. 1.5 Gridded Coverslip, 14 mm Glass Diameter |
| P35G-1.5-14-CGRD-D | No. 1.5 Gridded Coverslip Facing Down, 14 mm Glass Diameter |



50 mm MatTek Dishes

Low sidewalls make 50 mm dishes ideal for micro-injection. 50 mm dish covers snap securely to the bottom, making them perfect for atmosphere maintenance.

PART NO.

DESCRIPTION

50 mm Dish | No. 0 Coverslip

14 mm Glass Diameter

P50G-0-14-F

No. 0 Coverslip, 14 mm Glass Diameter

30 mm Glass Diameter

P50G-0-30-F

No. 0 Coverslip, 30 mm Glass Diameter

50 mm Dish | No. 1.5 Coverslip

14 mm Glass Diameter

P50G-1.5-14-F

No. 1.5 Coverslip, 14 mm Glass Diameter

P50GC-1.5-14-F

No. 1.5 Poly-D-Lysine Coated Coverslip, 14 mm Glass Diameter

30 mm Glass Diameter

P50G-1.5-30-F

No. 1.5 Coverslip, 30 mm Glass Diameter

50 mm Dish | No. 1.5 Gridded Coverslip

14 mm Glass Diameter

P50G-1.5-14-FGRD

No. 1.5 Gridded Coverslip, 14 mm Glass Diameter

60 mm MatTek Dishes



The 60 mm dishes provide an increased area for a greater volume of media useful for many cell culture applications, and also have low sidewalls for microinjection.

PART NO.

DESCRIPTION

60 mm Dish | No. 1.5 Coverslip

20 mm Glass Diameter

P60G-1.5-20-F

No. 1.5 Coverslip, 20 mm Glass Diameter

30 mm Glass Diameter

P60G-1.5-30-F

No. 1.5 Coverslip, 30 mm Glass Diameter



100 mm MatTek Dishes

The largest surface area available in a standard size dish, the 100 mm dishes support cell culture that require higher volumes of media.

PART NO.

DESCRIPTION

100 mm Dish | No. 1.5 Coverslip

30 mm Glass Diameter

P100G-1.5-30-F

No. 1.5 Coverslip, 100 mm Glass Diameter

ACCESSORIES

Coverslip Removal Fluid

50mL Bottle. For applications requiring removal of the coverslip from the dish, follow our protocol for easy removal of the glass coverslip from your MatTek dishes under Q5 on our FAQ page.

PART NO: P DCF OS 30





MatTek Glass Bottom Dish FAQ

HOW ARE GLASS BOTTOM DISHES TYPICALLY USED?

MatTek's glass bottom dishes are available uncoated or coated with poly-D-lysine or collagen. All dishes are gamma irradiated to ensure sterility. A general procedure for their use is as follows.

- Maintain sterility: Open dishes in a sterile environment (e.g. laminar flow hood).
- Pre-equilibrate dishes: Incubate the dishes with culture medium. Pipet 2-3 ml of medium into the 35 mm dishes or 3-4 ml into the 50 mm dishes and incubate at 37° C for 15 minutes.
- Add cell suspension to microwell: Remove the culture medium by aspiration and plate cells onto the glass surface. Pipet 250 µl of the cell suspension (cells suspended in culture medium) into the 10 mm diameter microwells, 500 µl of cell suspension into the 14 mm microwells, or 1 ml of cell suspension into the 20 mm wells. Incubate the dishes for 1 hour at 37° C.
- Add additional medium: After 1 hour, gently fill the remainder of the dish with medium. Add 2-3 ml to the 35 mm dishes or 3-4 ml for the 50 mm dishes.

Note: Contact us for more information on detailed protocols for 60 mm , 100 mm dishes, or other sizes.

WHAT TYPE OF GLASS BOTTOM DISH SHOULD I USE TO GROW MY CELLS?

Many transformed or cancerous cell lines will grow on uncoated dishes. It is hard to predict which type of glass bottom dishes (uncoated, poly-d-lysine coated, or collagen coated) will work best with your specific cell type. Poly-D-lysine coated dishes work well for neuronal culture and for many primary cells; other cells prefer a collagen coating. Additionally, many researchers purchase our uncoated dishes and apply their own specialized coating. We offer complimentary technical support and can help you choose the best dish for your needs.

CAN THE COVERSIP BE REMOVED FROM THE GLASS BOTTOM DISHES?

Yes, but for most applications, cells grown in the glass bottom dish can be viewed without removal of the coverslip using a variety of inverted and upright microscopes. We also provide coverslip removal fluid that works with our dishes.

CAN GLASS BOTTOM DISHES BE RE-USED?

MatTek Glass Bottom Dishes are meant for single-use experiments. We do NOT recommend re-using the glass bottom dishes. The re-use of dishes will introduce uncontrolled variables into your experiments which may affect the phenomenon under study. MatTek Dishes are also made of polystyrene and cannot be autoclaved.

Glass Bottom Multi-Well Culture Plates

MatTek's glass bottom multi-well plates combine the highest quality glass with the ability to grow up to 384 cultures under identical conditions.



For most applications, a 14 mm glass microwell is sufficient. Choose the 20 mm diameter microwell to maximize viewing area, or a 5 mm, 6 mm, or 10 mm diameter if your application is cell-, media-, or reagent-limited.

STERILE

All MatTek dishes are sterilized with gamma irradiation

MICROWELL APERTURE

| | 2.78 mm | 5 mm | 6 mm | 10 mm | 13 mm | 14 mm | 20 mm |
|----------|---------|------|------|-------|-------|-------|-------|
| 6-well | | | | ■ | | ■ | ■ |
| 12-well | | | | ■ | | ■ | |
| 24-well | | | | ■ | ■ | | |
| 48-well | | | ■ | | | | |
| 96-well | | ■ | | | | | |
| 384-well | ■ | | | | | | |

For most applications, No. 1.5 is the preferred coverslip thickness, especially for optimizing image quality with high numerical aperture objectives. The No. 0 coverslip gives you the most working distance and may be useful for thicker specimens.

GLASS THICKNESS

| | No. 0 | No. 1.0 | No. 1.5 |
|----------------|-------|---------|---------|
| 6-well | ■ | ■ | ■ |
| 12-well | ■ | ■ | ■ |
| 24-well | ■ | ■ | ■ |
| 48-well | | | ■ |
| 96-well | | | ■ |
| 96-well black | | | ■ |
| 384-well black | | | ■ |

Coatings: MatTek offers some poly-D-lysine and collagen-coated dishes as standard stocked items, but unique configurations may be custom-made as special order items.

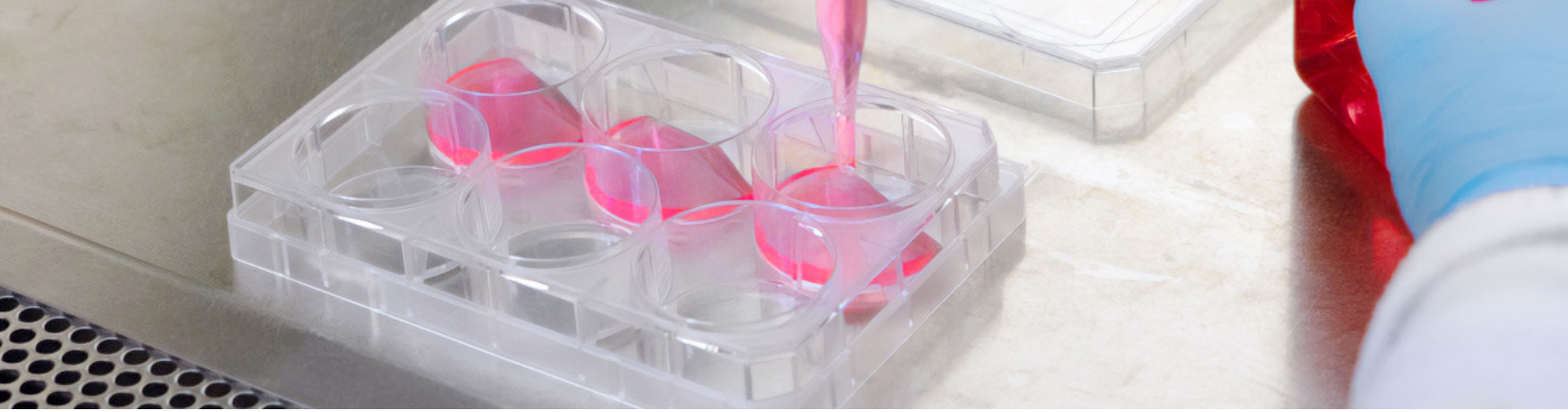


Culture up
to 384
samples in
our
glass
bottom
multi-well
plates

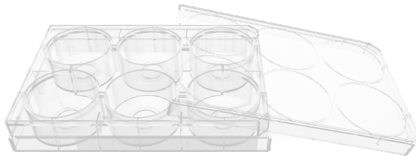
WHAT ARE THE ADVANTAGES OF USING GLASS BOTTOM MULTI-WELL PLATES COMPARED TO THE STANDARD GLASS BOTTOM CULTURE DISHES?

- The main advantage of the glass bottom multi-well plates is the ability to grow 6, 12, 24, 48, 96 or 384 cultures under identical conditions in the same culture plate. The glass bottom multi-well plates are ideal for high throughput and high-content screening applications.
- Analysis using the glass bottom multi-well plates is streamlined because only one plate (versus multiple Petri dishes) needs to be handled.
- For a number of applications, treatment of the cultures (e.g., irradiation) is simplified using glass bottom multi-well plates.
- Smaller wells in the glass bottom multi-well plates allow for decreased volumes of precious reagents.
- Compatibility with existing common laboratory equipment like multi-channel pipettes and plate readers.



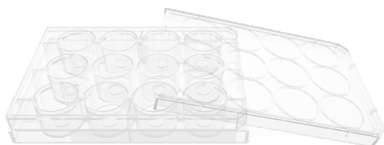


Glass Bottom 6-well Plates



Available in 10 mm, 14 mm, or 20 mm glass diameters, these 6-well plates offer the largest culture surface of any of our glass bottom multi-well plates.

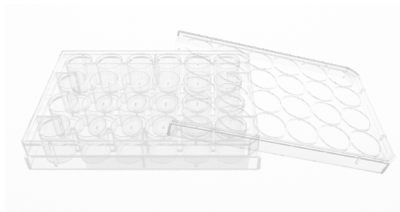
| PART NO. | DESCRIPTION |
|---|---|
| 6-well Plate No. 0 Coverslip | |
| 10 mm Glass Diameter | |
| P06G-0-10-F | No. 0 Coverslip, 10 mm Glass Diameter |
| 14 mm Glass Diameter | |
| P06G-0-14-F | No. 0 Coverslip, 14 mm Glass Diameter |
| 20 mm Glass Diameter | |
| P06G-0-20-F | No. 0 Coverslip, 20 mm Glass Diameter |
| 6-well Plate No. 1 Coverslip | |
| 10 mm Glass Diameter | |
| P06G-1.0-10-F | No. 1.0 Coverslip, 10 mm Glass Diameter |
| 14 mm Glass Diameter | |
| P06G-1.0-14-F | No. 1.0 Coverslip, 14 mm Glass Diameter |
| 20 mm Glass Diameter | |
| P06G-1.0-20-F | No. 1.0 Coverslip, 20 mm Glass Diameter |
| 6-well Plate No. 1.5 Coverslip | |
| 10 mm Glass Diameter | |
| P06G-1.5-10-F | No. 1.5 Coverslip, 10 mm Glass Diameter |
| 14 mm Glass Diameter | |
| P06G-1.5-14-F | No. 1.5 Coverslip, 14 mm Glass Diameter |
| 20 mm Glass Diameter | |
| P06G-1.5-20-F | No. 1.5 Coverslip, 20 mm Glass Diameter |



Glass Bottom 12-well Plates

Available in 10 mm or 14 mm glass diameters, these 12-well plates are perfect for testing in triplicate.

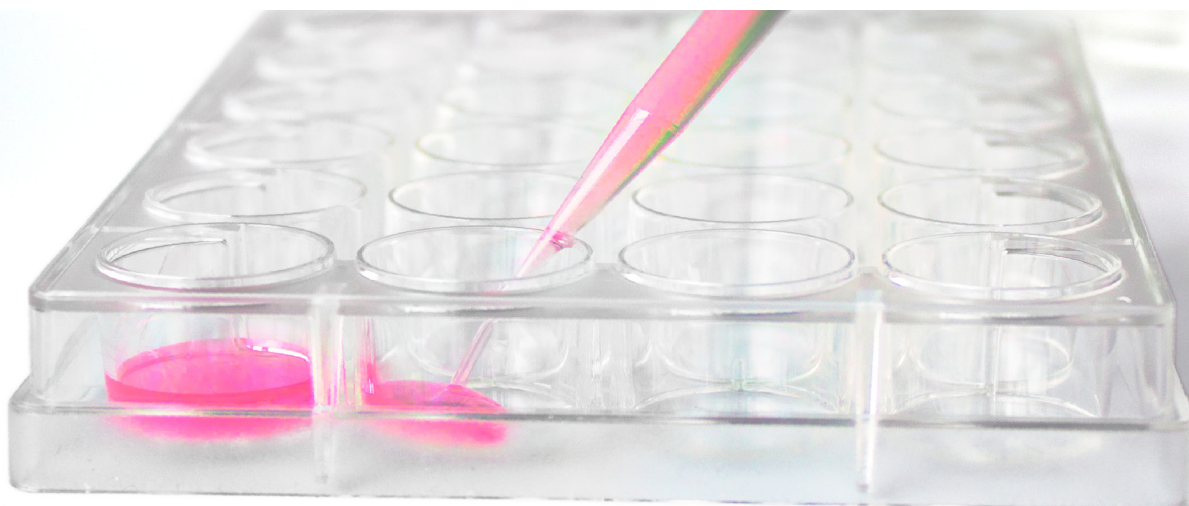
| PART NO. | DESCRIPTION |
|--|---|
| 12-well Plate No. 0 Coverslip | |
| 10 mm Glass Diameter | |
| P12G-0-10-F | No. 0 Coverslip, 10 mm Glass Diameter |
| 14 mm Glass Diameter | |
| P12G-0-14-F | No. 0 Coverslip, 14 mm Glass Diameter |
| 12-well Plate No. 1 Coverslip | |
| 10 mm Glass Diameter | |
| P12G-1.0-10-F | No. 1.0 Coverslip, 10 mm Glass Diameter |
| 14 mm Glass Diameter | |
| P12G-1.0-14-F | No. 1.0 Coverslip, 14 mm Glass Diameter |
| 12-well Plate No. 1.5 Coverslip | |
| 10 mm Glass Diameter | |
| P12G-1.5-10-F | No. 1.5 Coverslip, 10 mm Glass Diameter |
| 14 mm Glass Diameter | |
| P12G-1.5-14-F | No. 1.5 Coverslip, 14 mm Glass Diameter |

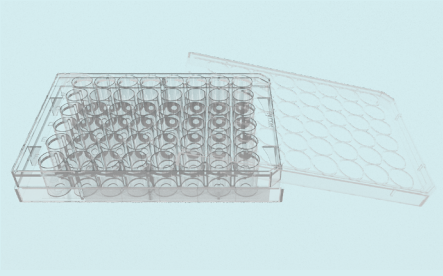


Glass Bottom 24-well Plates

Our 24-well plates are available in 10 mm or 13 mm microwell diameters.

| PART NO. | DESCRIPTION |
|--|---|
| 24-well Plate No. 0 Coverslip | |
| 10 mm Glass Diameter | |
| P24G-0-10-F | No. 0 Coverslip, 10 mm Glass Diameter |
| 13 mm Glass Diameter | |
| P24G-0-13-F | No. 0 Coverslip, 13 mm Glass Diameter |
| 24-well Plate No. 1.0 Coverslip | |
| 10 mm Glass Diameter | |
| P24G-1.0-10-F | No. 1.0 Coverslip, 10 mm Glass Diameter |
| 13 mm Glass Diameter | |
| P24G-1.0-13-F | No. 1.0 Coverslip, 13 mm Glass Diameter |
| 24-well Plate No. 1.5 Coverslip | |
| 10 mm Glass Diameter | |
| P24G-1.5-10-F | No. 1.5 Coverslip, 10 mm Glass Diameter |
| 13 mm Glass Diameter | |
| P24G-1.5-13-F | No. 1.5 Coverslip, 13 mm Glass Diameter |
| 24-well Plate No. 1.5 Coverslip | |
| Multi-well Plate Cover | |
| P24GTOP-1.5-F | No. 1.5 Coverslip |





Glass Bottom 48-well Plates

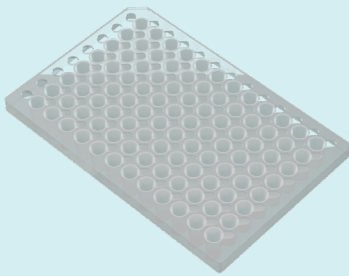
48-well plates are offered with a 6 mm glass microwell diameter.

| PART NO. | DESCRIPTION |
|----------|-------------|
|----------|-------------|

| | |
|--|--|
| 48-well Plate No. 1.5 Coverslip | |
|--|--|

6 mm Glass Diameter

| | |
|--------------|--|
| P48G-1.5-6-F | No. 1.5 Coverslip, 6 mm Glass Diameter |
|--------------|--|



Glass Bottom 96-well Plates

MatTek's signature high-quality cover glass is combined with the high-throughput convenience of a standard 96-well plate for brilliant imaging.

| PART NO. | DESCRIPTION |
|----------|-------------|
|----------|-------------|

| | |
|--|--|
| 96-well Plate No. 0 Coverslip | |
|--|--|

5 mm Glass Diameter

| | |
|------------|--------------------------------------|
| P96G-0-5-F | No. 0 Coverslip, 5 mm Glass Diameter |
|------------|--------------------------------------|

| | |
|--|--|
| 96-well Plate No. 1.5 Coverslip | |
|--|--|

5 mm Glass Diameter

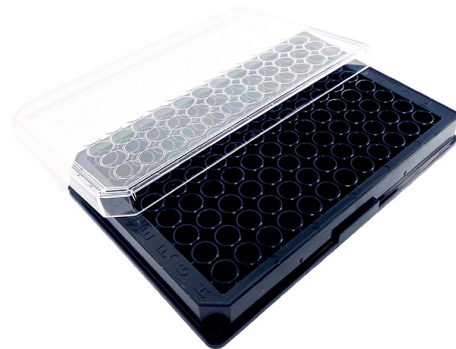
| | |
|----------------|--|
| P96G-1.5-5-F | No. 1.5 Coverslip, 5 mm Glass Diameter |
| P96GC-1.5-5-F | No. 1.5 Coverslip, 5 mm Glass Diameter, Poly-D-Lysine Coated |
| PBK96G-1.5-5-F | No. 1.5 Coverslip, 5 mm Glass Diameter, Black Plate |

FEATURED PRODUCT

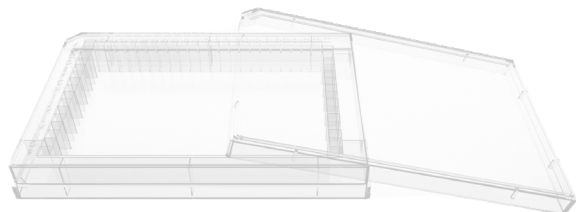
Glass Bottom 96-well Black Plates

Minimized backscattered light/background fluorescence. The glass bottom provides superior high-resolution imaging, while the black plate eliminates well-to-well crosstalk.

PART NO: PBK96G-1.5-5-F



Single Specimen Plate



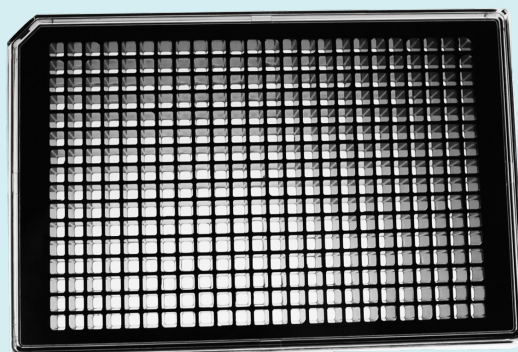
With a viewing area of 98mm X 67mm, MatTek's single specimen plate provides the greatest possible culture surface in a standard size plate. This generous viewing area makes it ideal for whole-animal or whole-organ imaging.

| PART NO. | DESCRIPTION |
|---|-------------|
| Single Specimen Plate No. 1.5 Coverslip Uncoated | |

| | |
|-------------------|---|
| P384G-1.5-10872-C | No. 1.5 Coverslip, 98mm x 67mm viewing area |
|-------------------|---|

Glass Bottom 384-well Black Plates

MatTek's glass bottom 384-well plates are ideal for high-throughput performance applications where low autofluorescence and exceptional optical clarity are needed. Ideal for high-resolution imaging, sensitive fluorescence, and confocal microscopy such as single molecule detection (SMD). Transmission measurements are capable in the wavelength range above 340nm.



| PART NO. | DESCRIPTION |
|--|-------------|
| 384-well Plate No. 1.5 Coverslip Uncoated | |

| | |
|---------------|--------------------------------|
| PBK384G-1.5-C | No. 1.5 Coverslip, Black Plate |
|---------------|--------------------------------|

Coverslips and Coverslip Kits

The highest quality German borosilicate glass coverslips in a wide range of sizes and thicknesses. Fix or stain your samples and finish with MatTek coverslips for immaculate slides and brilliant images or purchase our Coverslip Kits for the convenience of coverslips and Petri dishes in one package.



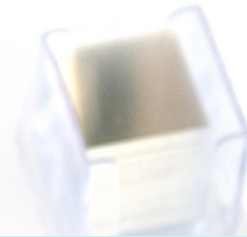
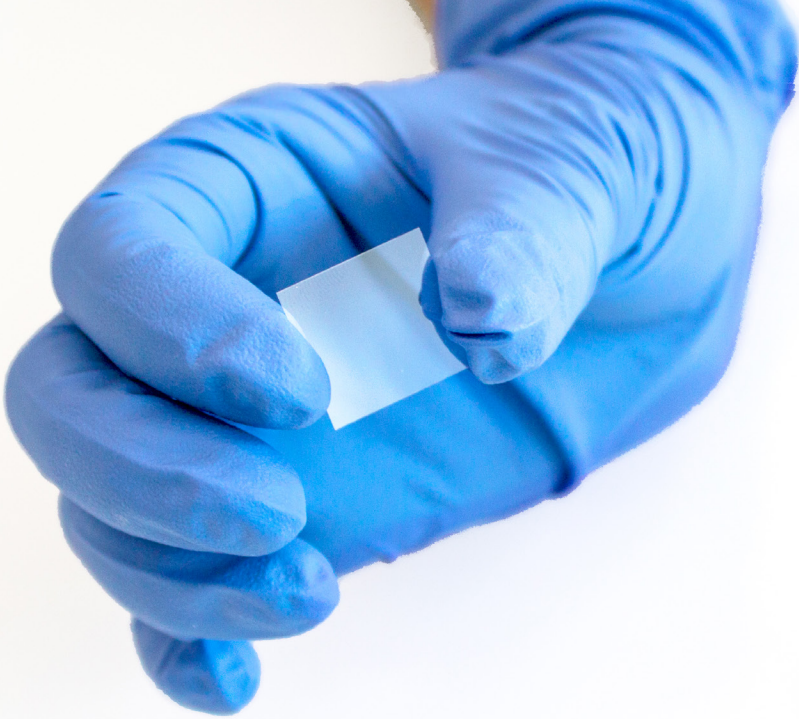
Select coverslips in square (gridded), rectangular, or round.

GLASS COVERSIP KITS

Ideal for amniocentesis, solid tumors, and chorionic villus samples (CVS), MatTek coverslip kits come pre-cleaned and sterilized. Our products are gamma-irradiated and come with a sterility guarantee.

GLASS COVERSGLIPS

The highest quality German borosilicate glass coverslips in a wide range of sizes and thicknesses.



Coverslip Kits

Ideal for amniocentesis, solid tumors, and chorionic villus samples (CVS), MatTek Coverslip Kits come pre-cleaned and sterilized. Our products are gamma-irradiated and come with a sterility guarantee.

PART NO.

DESCRIPTION

No. 1.5 Coverslip

35 mm Glass Diameter

| | |
|----------|--|
| CSGK/F | No. 1.5 Coverslip (22 x 22 mm), 35 mm Falcon® Dish |
| CSGK/F-F | Same as CSGK/F, also contains foam protection |
| CSGK/M | No. 1.5 Coverslip (22 x 22 mm), 35 mm MatTek Dish |
| CSGK/N | No. 1.5 Coverslip (22 x 22 mm), 35 mm Nunc® Dish |



Coverslips

High-quality glass in a wide range of sizes and thicknesses to suit your needs.

PART NO.

DESCRIPTION

Rectangular Coverslips

| | |
|---------------|---|
| PCS-1.5-5024 | 50 mm x 24 mm Coverslip, No. 1.5 Glass Thickness |
| PCS-0-10872 | 108 mm x 72 mm Coverslip, No. 0 Glass Thickness |
| PCS-1.5-10872 | 108 mm x 72 mm Coverslip, No. 1.5 Glass Thickness |

15 mm Round Coverslips

| | |
|------------|--|
| PCS-0-15 | 15 mm Coverslip, No. 0 Glass Thickness |
| PCS-1.0-15 | 15 mm Coverslip, No. 1.0 Glass Thickness |
| PCS-1.5-15 | 15 mm Coverslip, No. 1.5 Glass Thickness |

17 mm Round Coverslips

| | |
|------------|--|
| PCS-0-17 | 17 mm Coverslip, No. 0 Glass Thickness |
| PCS-1.0-17 | 17 mm Coverslip, No. 1.0 Glass Thickness |
| PCS-1.5-17 | 17 mm Coverslip, No. 1.5 Glass Thickness |

18 mm Round Coverslips

| | |
|------------|--|
| PCS-0-18 | 18 mm Coverslip, No. 0 Glass Thickness |
| PCS-1.0-18 | 18 mm Coverslip, No. 1.0 Glass Thickness |
| PCS-1.5-18 | 18 mm Coverslip, No. 1.5 Glass Thickness |

18 x 18 mm Square Coverslips

| | |
|------------------|--|
| PCS-0-1818 | 18 x 18 mm Coverslip, No. 0 Glass Thickness |
| PCS-1.0-1818 | 18 x 18 mm Coverslip, No. 1.0 Glass Thickness |
| PCS-1.5-1818 | 18 x 18 mm Coverslip, No. 1.5 Glass Thickness |
| PCS-170-1818 | 18 x 18 mm Coverslip, No. 1.5 (High Tolerance) Glass Thickness |
| PCS-1.5-1818-GRD | 18 x 18 mm Gridded Coverslip, No. 1.5 Glass Thickness |

22 x 22 mm Square Coverslips

| | |
|--------------|---|
| PCS-1.5-2222 | 22 x 22 mm Coverslip, No. 1.5 Glass Thickness |
|--------------|---|

25 mm Round Coverslips

| | |
|------------|--|
| PCS-0-25 | 25 mm Coverslip, No. 0 Glass Thickness |
| PCS-1.0-25 | 25 mm Coverslip, No. 1.0 Glass Thickness |
| PCS-1.5-25 | 25 mm Coverslip, No. 1.5 Glass Thickness |

35 mm Round Coverslips

| | |
|------------|--|
| PCS-0-35 | 35 mm Coverslip, No. 0 Glass Thickness |
| PCS-1.5-35 | 35 mm Coverslip, no. 1.5 Glass Thickness |

Chambered Cell Culture Slides

MatTek's Chambered Cell Culture Slides offer the ability to culture up to 8 different conditions on a single glass microscope slide. Culture, stain and examine your samples under a microscope without cell transfer.

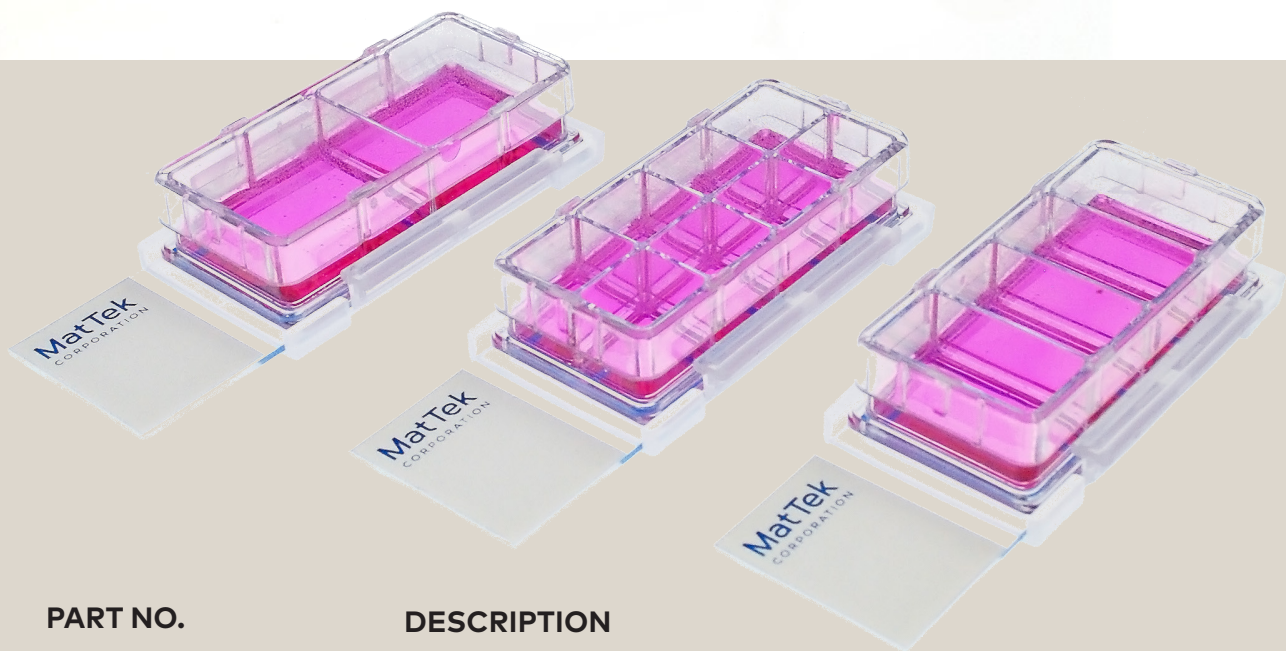


Unlike similar products, MatTek's Culture Slides do not use any adhesives, instead, they have a unique "snap-off" chamber apparatus that can be removed without the need for special tools. Naturally charged glass provides a superior growth surface for adherent cells.

STERILE

All MatTek Chambered Cell Culture Slides are sterilized with gamma irradiation





| PART NO. | DESCRIPTION |
|-----------------|--------------------|
|-----------------|--------------------|

| | |
|--|--|
| 2-well Chambered Cell Culture Slide | |
|--|--|

| | |
|-------|------------|
| CCS-2 | 2-well CCS |
|-------|------------|

| | |
|--|--|
| 4-well Chambered Cell Culture Slide | |
|--|--|

| | |
|-------|------------|
| CCS-4 | 4-well CCS |
|-------|------------|

| | |
|--|--|
| 8-well Chambered Cell Culture Slide | |
|--|--|

| | |
|-------|------------|
| CCS-8 | 8-well CCS |
|-------|------------|

PermaCell Cell Culture Inserts

MatTek PermaCell Inserts are advanced cell culture devices used for the growth and differentiation of cells. The 8mm inserts are uniquely made to function as either standing or hanging inserts, allowing researchers increased flexibility when designing studies. The 96-well insert plates consist of an array of 96 membrane coated wells in a single tray and allow for high-throughput processing.



The 8 mm inserts are sold as individual components in 24-well plates, and can be used in 6-, 12-, and 24-well culture plate formats. Various membrane types allow researchers to select inserts for specific needs, such as cell attachment, culture, differentiation, drug transport/permeability, and imaging studies. PermaCell Inserts can also be purchased as 96-well insert plates which consist of an array of 96 membrane wells connected into a single, rigid tray for easier handling and feeding

STERILE

All MatTek PermaCell Inserts are sterilized with gamma irradiation

MEMBRANE TYPES

- Polycarbonate (PC)
- Polyethylene terephthalate (PET)
- Teflon (PTFE)
- Specialty membranes also available

| Membrane Type | Pore Size | Pore Density | Transparent in Culture | Insert Size | Count | Part No. |
|---------------|-----------|---|------------------------|-------------|------------|----------------|
| PET | 0.4 µm | 1 x 10 ⁸ pores/cm ² | No | 24-well | 24 inserts | CCI24-PET-0.4 |
| PTFE | 0.4 µm | 1 x 10 ⁸ pores/cm ² | Yes | 24-well | 24 inserts | CCI24-PTFE-0.4 |
| PC | 0.4 µm | 1 x 10 ⁸ pores/cm ² | No | 24-well | 24 inserts | CCI24-PC-0.4 |
| | 8 µm | 1 x 10 ⁵ pores/cm ² | Yes | 24-well | 24 inserts | CCI24-PC-8 |

8 MM INSERT SPECIFICATIONS

| | |
|-------------------------|---|
| Height (including feet) | 8.8 mm |
| Height of Feet | 0.8 mm |
| Outer Diameter | 13.6 mm |
| Inner Diameter | 8.8 mm |
| Membrane Area | 0.6 cm ² |
| Membrane Pore Sizes | 0.4, 8 µm |
| Solvent Compatibility | Same as polystyrene |
| Sterility | Gamma Irradiation, shipped sterile 24-well plates |
| Membrane Types | Polytetrafluorethylene (PTFE), Polyethylene terephthalate (PET), Polycarbonate (PC) |

96-WELL INSERT PLATE SPECIFICATIONS

| | |
|-----------------------|------------------------------------|
| Dimensions | 96-well Insert Plate |
| Receiver Plates | 127.8 x 85.5 mm |
| Well Depth | 12 mm |
| Membrane Diameter | 4.29 mm |
| Membrane Area | 0.11 cm ² |
| Membrane Pore Size | 0.4 µm |
| Solvent Compatibility | Same as polystyrene |
| Sterility | Gamma Irradiation, shipped sterile |
| Membrane Types | Polyethylene terephthalate (PET) |



PermaCell Inserts

MatTek PermaCell Inserts have a pore size of 0.4 μm or 8 μm and are used for cell attachment, cell culture, cell differentiation, drug transport, and permeability studies.

These 8 mm inserts come packaged and pre-inserted in a 24-well plate. The 96-well insert plates come packaged and pre-inserted in a 96-well receiver plate.

| PART NO. | DESCRIPTION |
|----------|-------------|
|----------|-------------|

PermaCell Inserts | 0.4 μm

| | |
|-------------------|--|
| CCI24-PET-0.4 | 24-well, 0.4 μm , 24 Polyethelyene Terephthalate (PET) Membrane Inserts |
| CCI24-PTFE-0.4 | 24-well, 0.4 μm , 24 Polytetrafluorethylene (PTFE) Membrane Inserts |
| CCI24-PC-0.4 | 24-well, 0.4 μm , 24 Polycarbonate (PC) Membrane Inserts |
| CCI24-PTFE-CL-0.4 | 24-well, 0.4 μm , 24 Clear Polytetrafluorethylene (PTFE) Membrane Inserts |

PermaCell Inserts | 8 μm

| | |
|------------|---|
| CCI24-PC-8 | 24-well, 8 μm , 24 Polycarbonate (PC) Membrane Inserts |
|------------|---|

PermaCell Plates

MatTek PermaCell Insert Plates have a pore size of 0.4 μm and are used for high-throughput cell attachment, cell culture, cell differentiation, drug transport, and permeability studies. This 96-well insert plate design contains 96 inserts connected as a single, rigid tray and is compatible with TEER and robotics. The 96-well insert plates come packaged and pre-inserted in a 96-well receiver plate with lid. Single well reservoir plate may be purchased separately.



| PART NO. | DESCRIPTION |
|----------|-------------|
|----------|-------------|

96-well PermaCell Insert Plates | 0.4 μm


| | |
|---------------|--|
| CCI96-PET-0.4 | 96-well, 0.4 μm , 24 Polyethelyene Terephthalate (PET) Membrane |
|---------------|--|

96-well PermaCell Reciever Plates

| | |
|---------------|---|
| CCI96-RCVR-1 | Receiver plate with lid for the CCI96 plates with common (single) reservoir |
| CCI96-RCVR-96 | Receiver plate for CCI96 plates with 96 individual wells |

PermaCell Hang Top Lid

| | |
|---------------|--|
| CCI24-HANGTOP | 24-well plate with specialized lid for air-lifting of CCI24 PermaCell cell culture inserts (in a MatTek 24-well plate) |
| CCI12-HANGTOP | 12-well plate with specialized lid for air-lifting of CCI24 PermaCell cell culture inserts (in a MatTek 12-well plate) |

JOIN THE
BIO-CONVERGENCE
REVOLUTION 

MATTEK 

A BICO COMPANY

USA Headquarters

200 Homer Ave
Ashland, MA 01721
T: +1-508-881-6771

MatTek Europe
Mlynske Nivy 73
Bratislava 821 05 Slovak Republic
T: +421-2-3260-7401