

3D Microtissues, Primary Cells,
and Cultureware for Better Science

MATTEK

Now Part of Sartorius

We are Mattek, now a part of Sartorius.

Mattek Life Sciences was founded in 1985 and began producing human tissue models as reliable replacements for animal testing. Mattek's advanced three-dimensional (3D) tissue models of the skin, eye, oral, respiratory, vaginal, and intestinal systems empower academia and companies in the cosmetics, chemical, pharmaceutical, and consumer goods industries to achieve our goals of offering non-animal testing platforms that lower preclinical costs and provide more human-relevant results.

In addition to tissue models, Mattek offers a library of primary cells, customizable culture media, and a collection of cultureware products. Mattek's cultureware products are a valuable complement to our scientists. Available in standard dish and plate sizes, with the added feature of optically optimized surfaces suitable for high-quality imaging allows for experimental design convenience.

Mattek offers training for individuals interested in practicing various methods, consulting on specific problems, gaining valuable insights, or exploring the use of *in vitro* models for research. Contact us for more information.

Our Solutions

Supporting Your Research with
Advanced *In Vitro* Life Science
Products and Services



3D Microtissue Models

Mattek Tissue Models are produced from primary human cells and provide a microphysiological platform to model highly relevant and predictive human biology.

- EpiAirway
- EpiAlveolar
- EpiCorneal
- EpiDerm
- EpiDermFT
- EpiIntestinal
- EpiIntestinalFT
- EpiOcular
- EpiGingival
- EpiOral
- EpiNasal
- EpiVaginal
- EpiKidney
- EpiLiver
- HUREL Liver
- MelanoDerm
- Melanoma
- Psoriasis



Explore 3D Microtissue Models

Cultureware

Mattek Dishes have been the industry standard for cell culture and spectacular imaging for nearly 30 years. Experience the optics of glass in a petri dish.

- Glass Bottom Dishes
- Multi-well Plates
- Cover Slips
- Cover Slip Kits
- Chambered Cell Culture Slides
- PermaCell Culture Inserts



Explore Cultureware

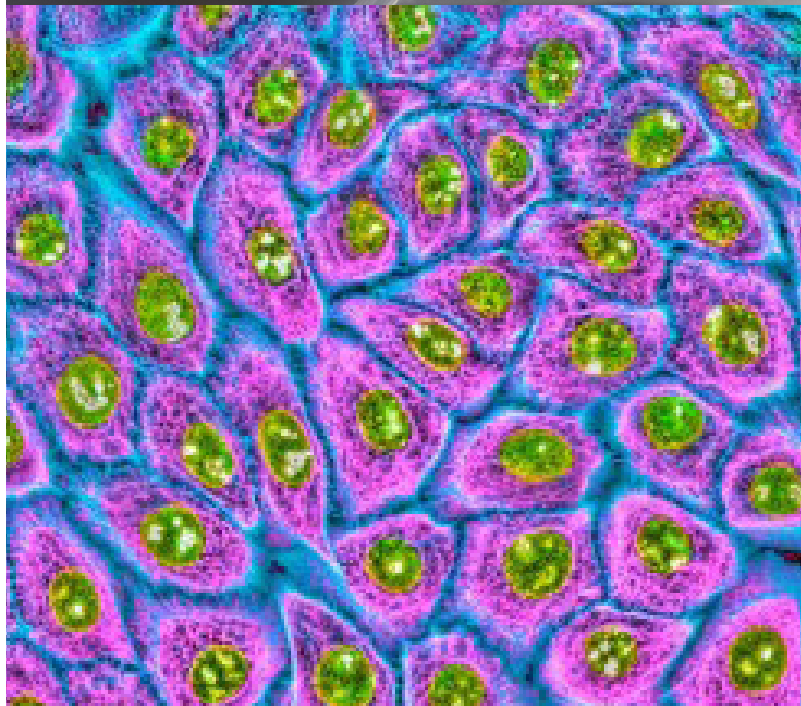
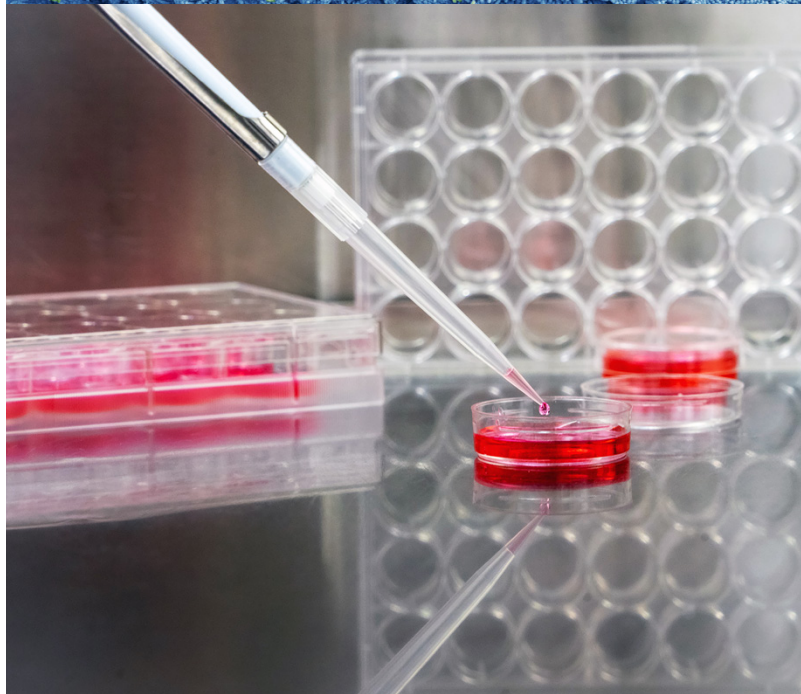
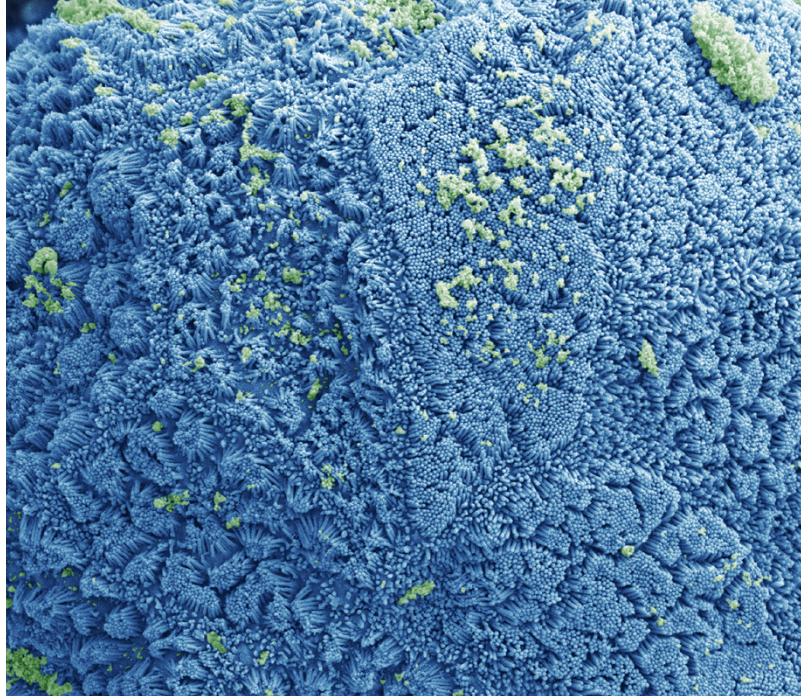
Primary Human Cells and Media

Mattek has been isolating cells for over 25 years, offering certified cells for 2D and 3D human biology.

- Skin Cells: Epidermal Keratinocytes, Melanocytes, Dermal Fibroblasts
- Airway Cells: Bronchial Epithelial Cells, Bronchial Fibroblasts
- Immune Cells: Dendritic Cells, Plasmacytoid Dendritic Cells, T-Cells

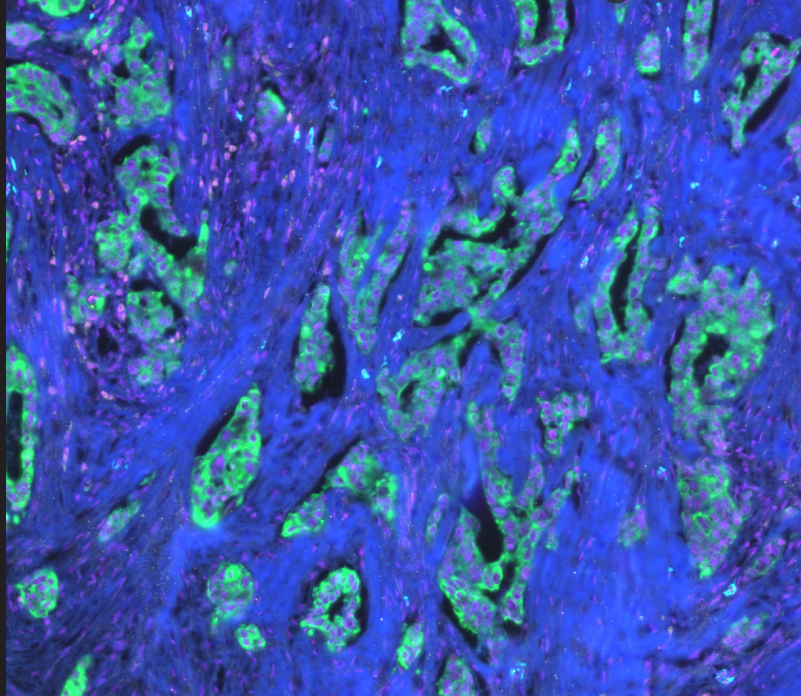


Explore Primary Cells and Media



Microtissue Toxicology Testing Services

Testing Services include advanced imaging and digital pathology tools for regulatory toxicology and efficacy studies across cosmetics, pharmaceutical, chemical, and household product industries. Their tissue models meet OECD and EU testing guidelines and are widely used in the US, Europe, and Asia.



VISIKOL® Contract Research and Testing Services:

Providing end-to-end services that include 3D tissue imaging, multiplex tissue imaging, digital pathology, high content imaging, 2D cell culture assays, 3D cell culture assays and *ex vivo* tissue slice assays. VISIKOL's expertise lies in both transforming tissues into actionable insights as well as bridging the gap between *in vitro* assays and *in vivo* results through the use of best-in-class cell culture models.

Liver diseases (e.g., NASH, NAFLD, HBV) are some of the most common diseases in the developed world. Liver toxicity is one of the most common sources of late-stage clinical failure.

▪ **In Vitro Liver, NASH, Fibrosis, and PCLS Services**

- HUREL® Micro Liver Model (2D Cell Culture)
- Primary Human Hepatocytes or HepaRG (Simple Spheroids)
- PHH or HepaRG with Non-Parenchymal Cells (Complex Spheroids)
- *Ex vivo* Tissue Slices

▪ **Blood Brain Barrier (BBB)**

- Immuno-oncology
- Histological Services
- Other tissue types:

- Kidney
- Lung

▪ **Other endpoints**

- Immunofluorescence
- 3D Tissue Imaging
- ELISA
- QPCR

▪ **More complex assays**

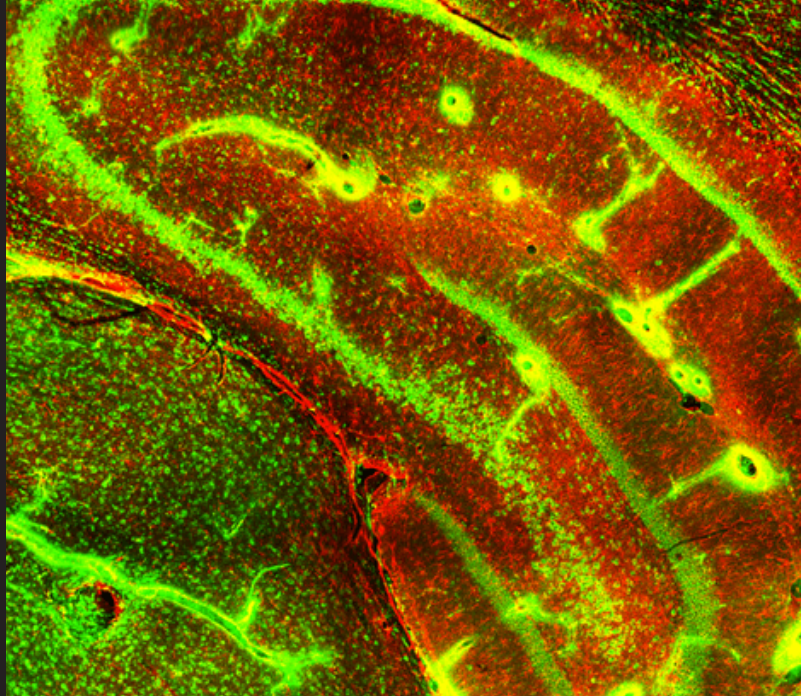
- Immune cell infiltration
- Therapeutic penetration
- Drug metabolism analysis (HPLC)

In addition to its services, VISIKOL® offers a suite of tissue clearing reagents and kits as well as the HUREL® Micro Liver portfolio of primary hepatocyte liver models.



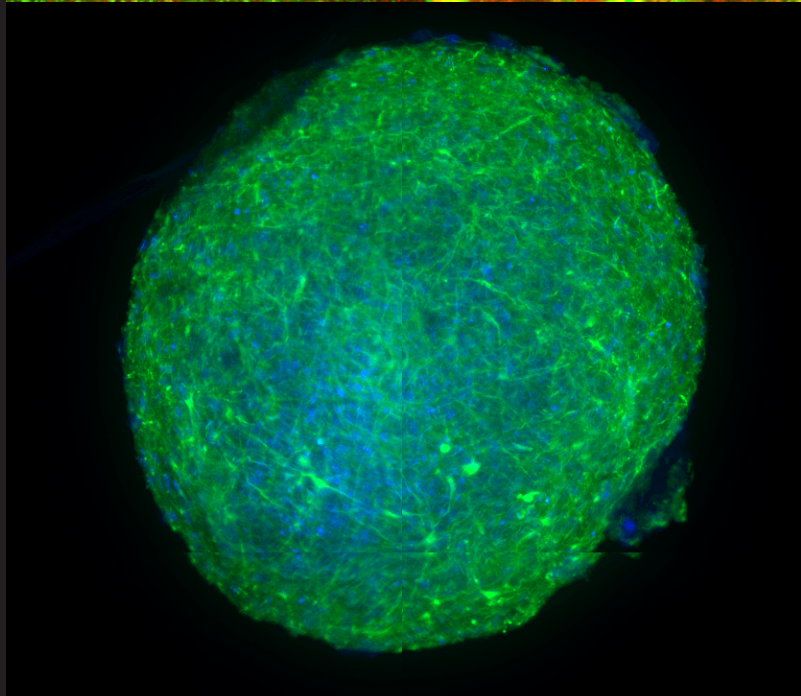
VISIKOL[®] HISTO[™]

A rapid, easy-to-use, affordable and reversible tissue clearing technique for whole tissues. Easily pair fluorescent protein and/or immunofluorescent labeling with tissue clearing for whole tissue (e.g. mouse brain, tumor) imaging using confocal, light sheet or 2-photon microscopy.



VISIKOL[®] HISTO-M[™]

A VISIKOL[®] HISTO[™] reagent designed specifically for use with 3D cell culture models (e.g. organoids, spheroids, microtissues) in high-throughput automated assays. The VISIKOL[®] HISTO-M[™] technology allows for complete 3D cell culture characterization instead of current assays that are biased towards only peripheral characterization.



VISIKOL[®] for Plant Biology[™]

A tissue clearing technique designed for use with plant samples for microscopic quality control. VISIKOL[®] for Plant Biology[™] is commonly used as a replacement for Hertwig's solution, which was previously used for this purpose and is a DEA controlled substance.



**Mattek North America
Headquarters**
200 Homer Ave
Ashland, MA 01721
US: + 1 (508) 881-6771

🌐 **For further information, visit**
www.mattek.com or email
support@mattek.com

Visikol, Inc.
200 Homer Ave
Ashland, MA 01721
US: + 1 (800) 615-8474

🌐 **For further information, visit**
www.visikol.com or email
info@visikol.com

**Mattek Europe
Mattek In Vitro Life Science
Laboratories**
Mlynské Nivy 73
821 05 Bratislava, Slovakia, EU
EU: +421 -2-3260-7401

🌐 **For further information, visit**
www.mattek.com or email
support@mattek.com

MATTEK

Now Part of Sartorius

VISIKOL

Now Part of Sartorius

MATTEK

Now Part of Sartorius

