

## Protocol

**TissueFab™ - GelMA-Conductive-Vis bioink, suitable for 3D bioprinting applications**

Protocol for Catalog No. [915963](#)

## Introduction

TissueFab™ - GelMA-Conductive-Vis bioink is a ready-to-use bioink which is formulated for electrical conductivity, high cell viability and printing fidelity. It is designed for extrusion-based 3D bioprinting and subsequent crosslinking with exposure to 405nm visible light. GelMA-Conductive bioinks can be used with most extrusion-based bioprinters, are biodegradable, and are compatible with human mesenchymal stem cells (hMSCs) and other diverse cell types. TissueFab™ - GelMA-Conductive-Vis bioink enables the precise fabrication of 3D cell models and tissue constructs for research in 3D cell biology, tissue engineering, in vitro tissue models, and regenerative medicine.

### Disclaimer

TissueFab™ - GelMA-Conductive-Vis bioink is for research use only; not suitable for human, animal, or other use. Please consult the Safety Data Sheet for information regarding hazards and safe handling practices.

## Specifications

<b>Storage</b>	Store TissueFab™ - GelMA-Conductive-Vis bioink at 2 - 8 °C. Protect from light by storing bottle in a foil bag or wrapping in aluminum foil.
<b>Stability</b>	Refer to the expiration date on the batch-specific Certificate of Analysis.

## Materials

### Materials supplied

The TissueFab™ - GelMA-Conductive-Vis bioink is supplied as follows:

Catalog Number	Quantity
<a href="#">915963</a>	1 × 10 mL bottle (1 unit)

### Materials required, but not supplied

- Cultured cells (visit our website for an up-to-date list of cell types) link: <https://www.sigmaaldrich.com/life-science/cell-culture/mammalian-cell-lines.html>
- Appropriate cell culture medium
- PBS (Cat. No. [D8537](#))
- Sterile pipette tips for transferring bioink
- Sterile printing cartridge, piston, and nozzle/needle for 3D printing
- Extrusion-based 3D bioprinter

- Water bath or incubator
- Micropipettes
- 405 nm light source

## Before you start: Important tips for optimal bioprinting results

---

**Optimize printing conditions.** Optimize printing conditions (e.g., nozzle diameter, printing speed, printing pressure, temperature, cell density) for the features of your 3D printer and for your application to ensure successful bioprinting. The suggestions below can guide you.

**Reduce bubble formation.** If the bioink has air bubbles, the bubbles may hamper bioprinting. Carefully handle the bioink when you mix and transfer it to avoid bubble formation. Do not vortex or shake vigorously.

**Aseptic techniques.** Follow standard aseptic handling techniques when you prepare and print the bioink, and during cell culture.

**Cell density.** Resuspend the cell pellet to the appropriate volume for the desired printed structure and cell density. Typical cell density for extrusion-based bioprinting is 1 to  $5 \times 10^6$  cells/mL. For example, Human bone marrow derived mesenchymal stem cells have been printed with TissueFab™ - GelMA-Conductive-Vis bioink at a concentration of  $5 \times 10^6$  cells/mL.

**Note:** The number of prints obtained from each 10-mL bottle of bioink (a unit) will vary depending on the structure that is printed. For example, each 10-mL bottle contains enough material to print a 30- $\mu$ L structure in each well of three 96-well plates or a 100- $\mu$ L structure in each well of four 24-well plates.

## Procedure

---

### A. Prepare bioink

1. Warm the 10-mL bottle of TissueFab™ - GelMA-Conductive-Vis bioink in a water bath or incubator set to 37 °C for 30 minutes or until the bioink becomes fluid, so that it is easy to pipette. The bioink is stable for at least 7 heating-cooling cycles.
2. When the bioink has become fluid, gently invert the TissueFab™ - GelMA-Conductive-Vis bioink bottle 10-15 times to make a homogeneous solution. DO NOT vortex or shake vigorously.

### B. Prepare bioink-cell solution

1. Centrifuge the cell suspension to obtain a cell pellet. Remove the supernatant carefully so that the cell pellet is not disrupted.
2. Resuspend the cell pellet at the desired cell density with the bioink solution by gently and slowly pipetting up and down several times. Ensure the cells are evenly distributed in the bioink solution by gently and slowly pipetting up and down several more times. Avoid creating air bubbles. DO NOT vortex or shake vigorously. Be careful not to dilute the bioink solution with cell culture medium because the medium might interfere with the printability of the bioink.
3. Pipette the bioink-cell solution into the desired printing cartridge. This step creates a filled printing cartridge.
4. Place the remaining bioink in a foil bag or wrap in aluminum foil and store at 4 °C to protect from heat and light.

### C. Bioprint

1. Cool the filled printing cartridge to 15-20 °C using a “temperature-controlled printhead”, if available, or place the cartridge in 4 °C refrigerator for 10–15 minutes to induce gelation. If your bioprinter has a temperature-controlled print bed, set temperature to 20 °C.
2. Follow the manufacturer’s 3D printer instructions. Load the print cartridge onto the 3D printer and print directly onto a Petri dish or into multi-well plates. Adjust the flow rate according the nozzle diameter, printing speed, printing pressure, and temperature.

#### **Example**

*Printer: Cellink BIO X™ or Cellink INKREDIBLE™ printer*

*Temperature: 20 °C*

*Flow rate (speed): 10 mm/s*

*Nozzle: 22G TT tapered needle*

*Pressure: 70-80 kPa*

### D. Crosslink

Place the light source directly above the 3D-bioprinted structure and expose the structure to 405 nm light (recommended settings: wavelength – 405 nm; power – 800 mW/cm<sup>2</sup>; distance – 8 cm; exposure – 30s). Use the appropriate distance and exposure time based on your light source. For 405 nm light sources usually available in desktop bioprinters, such as Cellink™ bioprinters (Bio X™ and INKREDIBLE™ printers,), distances of 3 cm or less and exposure times of 30-40 s or more are recommended.

The 3D-bioprinted structure is ready for culture or analysis immediately after crosslinking is done.

### E. Culture cells.

Culture the bioprinted tissue with the appropriate cell culture medium following standard tissue culture procedures.

## Troubleshooting

---

### **1. Bioink is incubated at 37°C for 30 minutes, but it is still gel.**

Possible reasons – Malfunction of incubator; bioink is crosslinked due to light exposure.

Solution – Make sure the temperature of incubator / water bath is correct and make sure the bioink bottle is properly and evenly heated in the incubator/water bath. Do not expose the bioink to light before printing.

### **2. Air bubble is trapped in the middle of bioink in the cartridge.**

Possible reason – Air bubble was created when the bioink was transferred (and/or was mixed with cells).

Solution - Warm the cartridge at 37°C for 5–10 minutes or until the bioink becomes fluid. Turn the cartridge so that the tip faces up to allow any air bubbles to exit from the tip of the cartridge. Gently tap the cartridge to help the air bubbles pass through the tip.

### 3. Printed structure spreads and does not hold its shape.

Possible reasons – Bioink was diluted with cell culture medium that remained in the cell pellet; bioink was not cooled sufficiently before printing; or the printing pressure is too high.

Solution – Do not dilute the bioink. Make sure the bioink has been cooled according to the instructions before printing. Adjust printing pressure to achieve sufficient flow of bioink.

### 4. Interrupted flow or no flow during printing.

Possible reason – Insufficient printing pressure or nozzle is partially or fully clogged.

Solution – Adjust the printing pressure to achieve sufficient flow of bioink. If the problem persists, change the nozzle.

### 5. Printed structure dissolves in cell culture medium.

Possible reason – Insufficient crosslinking; exposure to incorrect wavelength; malfunction of light source apparatus.

Solution – Make sure that the light source has proper wavelength and sufficient power output and that the printed structure is exposed to light according to the instructions.

## Related Products

Name	Cat. No.
TissueFab™ - GelAlg-Vis bioink	<a href="#">906913</a>
TissueFab™ - GelMA-UV bioink	<a href="#">905429</a>
TissueFab™ - GelMA-Vis bioink	<a href="#">906891</a>
TissueFab™ - Sacrificial bioink	<a href="#">906905</a>
TissueFab™ - GelMA-Conductive-UV bioink	<a href="#">915726</a>
TissueFab™ - GelMA-Bone-UV bioink	<a href="#">915025</a>
TissueFab™ - GelMA-Bone-Vis bioink	<a href="#">915033</a>

Copyright © 2021 Merck KGaA, Darmstadt, Germany and/or its affiliates. All rights reserved. Merck, the vibrant M, SigmaAldrich, and TissueFab are trademarks of Merck KGaA, Darmstadt, Germany or its affiliates. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly accessible resources. More information on our branded products and services on [MerckMillipore.com](https://www.MerckMillipore.com)

The life science business of Merck  
operates as MilliporeSigma in the  
U.S. and Canada.

**MERCK**