

## Product Information

### TrueGel3D Hydrogel Kits

CD cell-degradable crosslinker and RGD peptide

Catalog Number **TRUE1**

Storage Temperature  $-70\text{ }^{\circ}\text{C}$

## TECHNICAL BULLETIN

### Product Description

TrueGel3D hydrogel pre-configured kit is used to prepare a chemically defined hydrogel with stiffness of  $\sim 400\text{ Pa}$  (shear modulus) that can be customized to match that of native cell environment.

The TrueGel3D hydrogel is formed by crosslinking of RGD degradable polymer with CD cell-degradable crosslinker. The cells are encapsulated during crosslinking, where they can adhere to the polymer through the RGD peptide and grow within the hydrogel. The CD cell-degradable crosslinker is composed of matrix metalloprotease (MMP)-cleavable peptide (Pro-Leu-Gly-Leu-Trp-Ala), which allows cells to spread and migrate by secreting matrix metalloproteases (MMP1, MMP3, MMP7, and MMP9).

TrueGel3D hydrogel with RGD degradable polymer can be dissolved by treatment of TrueGel3D Enzymatic Cell Recovery Solution (Catalog Number TRUEENZ) to recover cells for post culture analysis.

### Components

RGD degradable polymer, lyophilized Catalog Number TRU-RGD	$2 \times 870\text{ }\mu\text{L}$
CD cell-degradable crosslinker, lyophilized Catalog Number TRU-CD	$200\text{ }\mu\text{L}$
TrueGel3D reconstitution Buffer Catalog Number TRU-REC	$2 \times 900\text{ }\mu\text{L}$
Water Catalog Number TRUWA	$600\text{ }\mu\text{L}$

### Precautions and Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Safety Data Sheet for information regarding hazards and safe handling practices.

### Preparation Instructions

RGD degradable polymer

Note: When stored at  $-70\text{ }^{\circ}\text{C}$ , warm RGD degradable polymer to room temperature before each use.

- Reconstitute a tube (RGD degradable polymer) with  $860\text{ }\mu\text{L}$  of buffer.
- Vortex until all material is dissolved.
- Incubate reconstituted polymer for 1 hour at room temperature.
- Centrifuge briefly; RGD degradable polymer is ready to use.

CD cell-degradable crosslinker

- Reconstitute a tube with  $188\text{ }\mu\text{L}$  of water to obtain concentration of  $20\text{ mM}$  thiol groups.
- Vortex until all material is dissolved.
- Incubate at room temperature for 5 minutes.
- Vortex and centrifuge the tube.
- CD cell-degradable crosslinker is ready to use.

### Storage/Stability

- The lyophilized powders may be stored unopened in the original bottles at  $-70\text{ }^{\circ}\text{C}$  for up to one year.
- Do not expose the crosslinker to air longer than necessary to avoid oxidation of thiol groups. After reconstitution, it can be stored at  $-20\text{ }^{\circ}\text{C}$  or  $-80\text{ }^{\circ}\text{C}$ .
- For short term storage ( $<1\text{ month}$ ) both RGD degradable polymer and reconstitution buffer are stored at  $4\text{ }^{\circ}\text{C}$ .
- Water can be stored between  $-70\text{ }^{\circ}\text{C}$  and room temperature.

## Procedures

### Formation of Hydrogel

All steps are performed in sterile hood.

1. Prepare cell suspension using culture medium, PBS, or any other physiological solution.  
Note: Cell suspension should be 20% of final volume of the gel.
2. Add RGD degradable polymer to cell suspension in a reaction tube and mix gently. Add CD cell-degradable crosslinker to the mixture and mix by pipetting few times. The volume ratio of each component is added as indicated in Table 1.

**Table 1.**

Gel Component Volumes

Components	Volume parts
RGD degradable polymer	18
Cell suspension	5
CD cell-degradable crosslinker	2
<b>Total gel volume</b>	<b>25</b>

3. Plate the mixture in a culture dish before it begins to solidify.  
Note: The mix will stay liquid between 1–4 minutes and afterward will begin to form a gel. Make sure to dispense all of the mix during the period after mixing
4. Incubate the mixture at 37 °C for 20 minutes. Alternatively, it can also be incubated at room temperature with a slightly longer incubation period.
5. Test gel formation by gently touching gel with pipette tip. It should not pull out threads of gels when retracting from the gel surface.
6. Add the culture medium to cover the gel.
7. Place the lid on culture dish and incubate for cell cultivation.
8. Replace the medium after 1 hour.
9. Change the medium as needed for proper growth of cells.

### Recovery of cells

TrueGel3D Enzymatic Cell Recovery Solution is used to dissolve the hydrogel matrix.

1. Add 300 µL of 1:20 diluted TrueGel3D Enzymatic Cell Recovery Solution to dissolve 25 µL of gel.  
Note: Rate of dissolution is increased if gels are cut into pieces
2. Incubate at 37 °C for 30–60 minutes.
3. Centrifuge the cell suspension and resuspend the pelleted cells in fresh medium or buffer.
4. Repeat step 3 twice to wash the remains of TrueGel3D Enzymatic Cell Recovery Solution from the gel components.
5. Cells are now ready to use for post culture analysis or to set up new hydrogel.  
Note: If TrueGel3D Enzymatic Cell Recovery Solution is not removed completely, it destabilizes the newly set up hydrogel.

### Reference

1. Knight, C.G. et al., A novel coumarin-labelled peptide for sensitive continuous assays of the matrix metalloproteinases. *FEBS Lett.*, **296**, 263–266 (1992).

BG,MAM 12/17-1