

Product Information

TrueGel3D Hydrogel Kits

SLO-PVA, PEG non cell-degradable crosslinker

Catalog Number **TRUE9**

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

TECHNICAL BULLETIN

Product Description

TrueGel3D Hydrogel Kits with SLO-PVA polymer are used to set up chemically defined, slow gelling hydrogels. It has higher stiffness that can be customized to match that of the native cell environment. The gel is formed by crosslinking of SLO-PVA polymers with PEG non cell-degradable crosslinkers. SLO-PVA polymers allow enough time to conveniently manipulate the solution unlike fast gelling hydrogels. Slow gelling hydrogels can be used in microchannels or syringes.

TrueGel3D Hydrogel with SLO-PVA polymer can be customized by adding TrueGel3D RGD peptide (Catalog Number TRUERGD) to provide attachment sites for cells. The cells are encapsulated during crosslinking, where they can adhere to the polymer through the RGD peptide and grow within the hydrogel.

Extracellular matrix (ECM) proteins (Fibronectin, Laminin) or other bioactive components like growth factors can also be added in the hydrogel mix: please refer to TrueGel3D Slow protocol online for more details.

Components

- SLO-PVA solution 170 μL
in phosphate buffer
Each tube contains 30 mmol/L reactive groups
Catalog Number TRU-SPVA
- PEG non cell-degradable crosslinker 200 μL
lyophilized
Each tube contains 20 mmol/L reactive groups
Catalog Number TRU-PEG
- TrueGel3D buffer 200 μL
10 \times concentrated, pH 7.2
Catalog Number TRU-B72
- Water 2 \times 1,500 μL
Catalog Number TRUWA

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

PEG non cell-degradable crosslinker

- Centrifuge the vial to make sure entire material is at the bottom of the tube.
- Add 188 μL of water to make a concentration of 20 mmol/L thiol groups.
- Vortex until all material is dissolved.
- Incubate at room temperature for 5 minutes
- Vortex and centrifuge the tube.
- PEG non cell-degradable crosslinker is now 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.
- SLO-PVA may be stored at $-70\text{ }^{\circ}\text{C}$ for long term and $4\text{ }^{\circ}\text{C}$ for short term.
- Do not expose the PEG non cell-degradable crosslinker/RGD peptide to air longer than necessary to avoid oxidation of thiol groups. After reconstitution, it can be stored at $-20\text{ }^{\circ}\text{C}$ or $-70\text{ }^{\circ}\text{C}$.
- Buffers are stored at $4\text{ }^{\circ}\text{C}$ for short term (<2 months) and for long term between $-20\text{ }^{\circ}\text{C}$ and $-70\text{ }^{\circ}\text{C}$.
- Water can be stored between $-70\text{ }^{\circ}\text{C}$ and room temperature.

Procedure**Formation of Hydrogel**

All steps are performed in sterile hood and the volume ratio of each component is added as indicated below.

| Components | Without peptide (µL) | With Peptide (µL) |
|---|----------------------|-------------------|
| Water | 16.6 | 15.3 |
| TrueGel3D buffer, 10× concentrated, pH 7.2 | 2.4 | 2.4 |
| SLO-PVA (30 mmol/L) | 2.0 | 2.5 |
| RGD peptide (20 mmol/L) | - | 0.8 |
| Cell suspension | 6.0 | 6.0 |
| PEG non cell-degradable crosslinker (20 mmol/L) | 3.0 | 3.0 |
| Total | 30.0 | 30.0 |

1. Prepare cell suspension using culture medium, PBS, or any other physiological solution.
2. Mix water, 10× TrueGel3D buffer, pH 7.2, and SLO-PVA in a reaction tube and mix well.
3. Add the RGD peptide (if applicable) to the reaction tube containing SLO-PVA and mix immediately to ensure homogenous distribution. Incubate for 20 min to allow attachment of the RGD peptide to the SLO-PVA polymer.
Note: If RGD peptide is not used, skip this step.
4. Add cell suspension to the reaction tube containing the polymer (SLO-PVA) to prepare cell suspension mix.
5. Add the PEG non cell-degradable crosslinker to the cell suspension mix and pipette a few times.
6. Incubate the solution for 10 minutes at room temperature.
Note: Do not incubate longer than 10 minutes as the solution will solidify and cannot be transferred through the pipette.
7. Resuspend cells by pipetting for a few times to ensure uniform distribution in the gel and transfer the solution in a sterile culture dish.
8. Incubate for 50 minutes at room temperature or at 37 °C.
9. Once gel has formed, add the cell culture medium until the gel is covered.
10. Incubate the culture dish in the incubator.
11. Replace the medium after 1 hour.
12. Change culture medium as required for proper growth of cells.

BG,MAM 01/18-1