

Product Information

TrueGel3D Enzymatic Cell Recovery Solution

Catalog Number **TRUEENZ**
Storage Temperature $-70\text{ }^{\circ}\text{C}$

TECHNICAL BULLETIN

Product Description

TrueGel3D Enzymatic Cell Recovery Solution (TRUEENZ) is used to dissolve the hydrogel matrix and recover cells for post culture analysis from dextran based hydrogels.

Dextranase isolated from *Chaetomium gracile* is the main constituent of TrueGel3D enzymatic cell recovery solution. It cleaves glycosidic bonds of dextran to recover both live and fixed cells without any cellular damage.

Component

TrueGel3D Enzymatic Cell Recovery Solution 500 μL

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.

Storage/Stability

TrueGel3D Enzymatic Cell Recovery Solution may be stored at $4\text{ }^{\circ}\text{C}$ for short term (<6 months), and at $-70\text{ }^{\circ}\text{C}$ for long term storage. Avoid frequent freeze-thaw cycles

Procedure

1. Add 300 μL of TrueGel3D Enzymatic Cell Recovery Solution diluted 1:20 with cell culture supernatant to dissolve 25 μL of gel.
Note: Increase rate of dissolution by cutting gels into small pieces.
2. Incubate at $37\text{ }^{\circ}\text{C}$ 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 remove the remaining TrueGel3D Enzymatic Cell Recovery Solution in 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 will destabilize the newly set up hydrogel

BG,MAM 12/17-1