



# Reproducible animal-free 3D cell culture.

## GrowDex hydrogels for reproducible 3D cell culture

GrowDex® hydrogels mimic the extracellular matrix and support cell growth and differentiation enabling convenient, reproducible and scalable animal-free culture of 3D spheroids and organoids. Available from Revvity, GrowDex hydrogels are composed of natural nanofibers of cellulose, extracted from wood sourced from sustainable and responsibly managed forests, and are a proven solution for automated cell culture and high-content screening.

### GrowDex®



GrowDex hydrogels

For research use only. Not for use in diagnostic procedures. GrowDex® is a registered trademark of UPM-Kymmene Corp.

#### Reproducible

Manufactured to the highest standards with strict quality control criteria to ensure lot to lot reproducibility.

#### Animal-free

No animal biomolecules interfering with readouts.

#### Convenient

Used and stored at room temperature, suitable for liquid handling systems and supplied in prepacked syringes.

Simple protocol - mix GrowDex hydrogels with media and cells.

#### Tunable

Optimal gel stiffness is achieved by dilution with cell culture media. Matrix composition can be optimized easily by supplementing cell culture media with specific biomolecules such as growth factors.

#### Optimized for imaging

Transparent, no auto-fluorescence, and tried and tested on Revvity high-content imaging systems.



## Product list

Description	Part number	Pack size
<b>GrowDex</b> The original hydrogel, highly biocompatible GrowDex has been used widely for culture of primary cells through to iPS and ES cells.	100103002	2.5 mL
	100103005	5 mL
	100103010	10 mL
<b>GrowDex-T</b> GrowDex-T brings superior imaging quality for phase contrast and brightfield imaging. It is completely transparent - ideal for imaging and automated analysis.	200103002	2.5 mL
	200103005	5 mL
	200103010	10 mL
<b>GrowDex-A</b> GrowDex -A is ideal if you need to customize your hydrogel. It has been modified by the addition of an avidin group so you can attach any biotinylated molecule.	300103002	2.5 mL
	300103005	5 mL
	300103010	10 mL
<b>GrowDase</b> GrowDase enzyme breaks down GrowDex hydrogels to form a solution in an easy, one-step, efficient cell recovery process.	900102001	1.25 mL
	900102002	2.5 mL

revvity