

Which GrowDex[®] hydrogel to choose?

GrowDex[®]

- For spheroid and organoid formation.
- **Examples: Cancer cells, lung, liver, stem cells**



Do you want to recover your cells from the matrix?
E.g. for downstream processes?

Yes, choose **GrowDase**



GrowDex[®]-T



Yes, choose **GrowDex-T**

Are you looking for exceptionally good imaging properties?

Yes

Do you also want to link molecules such as ECM proteins, growth factors or antibodies to the matrix?

GrowDex[®]-A



Yes, choose **GrowDex-A**

- For organoids and cells that require anchorage points, such as ECM-molecules.
- **Examples: Cancer cells, heart, muscle, intestine, vasculature**

- For spheroids and cells that have outgrowth or elongated morphology.
- **Examples: Neuronal cells, fibroblasts, MSCs**

FEATURE	GrowDex [®]	GrowDex [®] -T	GrowDex [®] -A
Cellulose type	Native	Anionic	Anionic
Cellulose fiber charge	Neutral	Negative	Negative
Stock concentration	1.5%	1.0%	1.0%
Working concentration range	0.2% - 1.5%	0.1% - 1.0%	0.1% - 1.0%
Visual appearance	Opaque	Transparent	Transparent
Brightfield	**	***	***
Phase contrast	**	***	***
Fluorecence	***	***	***
Ready-to-use	Yes	Yes	Yes, mix with biotinylated molecule
Shelf life	12 months	12 months	6 months

To order products online: www.revivity.com

To request a quote or place an order: www.revivity.com/contact-us

For technical support: LS.ReagentsTechSupport@revivity.com



Revivity, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (+1) 781 663 6900
www.revivity.com