

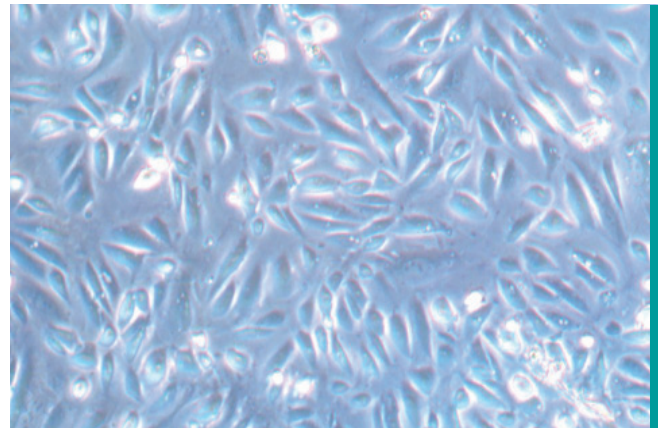
## Dissociated Tumor Cells (DTCs)

### What are DTCs?

Dissociated Tumor Cells (DTCs) are a single cell suspension that is dissociated from a solid tumor using enzymes and mechanical digestion. They have a heterogeneous cell composition.

Primary cells offer the closest known comparison to the original tissue and are key to research that focuses on advancing personalized medicine and drug development.

- Strict quality control to ensure the product meets specifications
- Complete clinical data set provided for each lot
- Cryopreserved to provide greater convenience and allow you to dictate your study schedule
- Sourced from our expansive ASTERAND™ Human Tissue collection network
- Availability of matched normal adjacent tissue and cells available



Renal Cell Carcinoma DTCs at 100x magnification.

## Personalized Medicine Applications

DTCs provide the closest representation of the clinical scenario. While no two patients' tumors are identical, there is a greater likelihood of a therapy being effective if it has been previously tested on DTCs from patients with similar health issues.

Researchers can use patient-derived DTCs to determine how specific cohorts of cancer cells will react to different drug therapies.

DTCs can also be incorporated into 3D micro-tumor arrays to provide highly efficient, high-throughput drug screening and chemosensitivity testing.

# Expertise In the Isolation and Propagation of Human Primary Cells

## Quality Control

Providing high-quality products is our top priority, a pathologist will review each tumor sample at collection to confirm that it is the correct tissue and organ site. Also, they will confirm the diagnosis matches the clinical data.

Our quality control processes ensure that:

- Lots are sterile and pathogen-free
- Cells grown to confluency in cell culture to validate that they are viable and will grow as expected



All information is detailed on the Certificate of Analysis accompanying the cell lot.

CANCER COHORTS	PRODUCT NUMBER	AVAILABLE DTC SIZES
Human Breast DTC	HPCDTCBR	<1 Million Cells/Vial
Human Kidney DTC	HPCDTCKID	1-2 Million Cells/Vial
Human Liver DTC	HPCDTCLIV	2-3 Million Cells/Vial
Human Lung DTC	HPCDTCLUN	3-4 Million Cells/Vial
Human Colon DTC	HPCDTCCRC	4-5 Million Cells/Vial
Human Ovarian DTC	HPCDTCOVA	>5 Million Cells/Vial

And more...



## Related Products:

- Primary Human Cancer Associated Fibroblasts
- Primary Human Cancer Associated Epithelial Cells
- SUM Breast Cancer Cell Lines

Visit [BioIVT.com/inventory](http://BioIVT.com/inventory) to view our ready-to-ship inventory.



### For Inquiries:

North America & Asia Pacific T +1 516 483 1196 F +1 516 483 4683 E [customerservice@bioivt.com](mailto:customerservice@bioivt.com)

Europe, Middle East & Africa T +44 (0) 1444 250010 F +44 (0) 1444 250066 E [cs europe@bioivt.com](mailto:cs europe@bioivt.com)