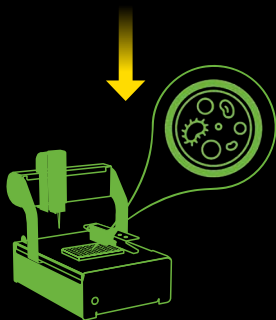


# FROM BIOLOGICS TO BURGERS: The Journey of Cultured Meat

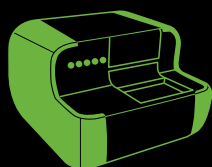
As the demand for sustainable and ethical meat sources grows, cultured meat offers a promising solution. Discover the fascinating journey of lab-grown meat from its inception in a petri dish to its culmination on our dinner plates. Lab-grown meat represents a blend of cutting-edge technology and our quest for sustainability. It promises to revolutionize our dinner plates, making them more ethical and eco-friendly without compromising the flavors we cherish.



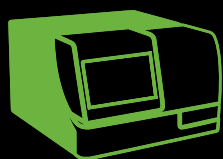
**Animal tissue collection:** The process starts with a small biopsy from a live animal, typically taken from muscle tissue. This is cruelty-free and doesn't harm the animal.



**Single-cell isolation:** Using the DispenceCell™ Single-Cell Dispenser, individual cells are isolated from the tissue. These cells have the potential to grow and differentiate into muscle tissues.



**Monoclonality verification and growth optimization:** The isolated cells are cultivated in a controlled environment. With the CloneSelect® Imager, the best single-cell clones are identified and further cultivated. This step ensures that the cells have the optimal growth characteristics.



**Characterization and safety assessment:** As the cells multiply and form muscle tissue, their growth and characteristics are monitored using the SpectraMax® iD5 Multi-Mode Microplate Reader. This ensures that the meat produced meets the desired quality and nutritional standards.



The magic of cultured meat transforms simple cell samples into flavorful feasts. Beginning with a tiny tissue biopsy, science and technology converge to cultivate cells, optimize their growth, and monitor their quality. The end result? A sustainable and ethical dish that satisfies both the palate and the planet.

For more information, visit  
[www.moleculardevices.com/cultured-meats](http://www.moleculardevices.com/cultured-meats)



## Contact Us

Phone: +1.800.635.5577  
Web: [www.moleculardevices.com](http://www.moleculardevices.com)  
Email: [Info@moldev.com](mailto:Info@moldev.com)  
Check our website for a current listing of worldwide distributors.

## Regional Offices

USA and Canada	+1.800.635.5577	Taiwan/Hong Kong	+886.2.2656.7585
United Kingdom	+44.118.944.8000	Japan	+81.3.6362.9109
Europe*	00800.665.32860	South Korea	+82.2.3471.9531
China	+86.4008203586	India	+91.73.8661.1198

\*Austria, Belgium, Denmark, Finland, France, Germany, Iceland, Ireland, Italy, Luxembourg, Netherlands, Portugal Spain, Sweden, Switzerland and United Kingdom