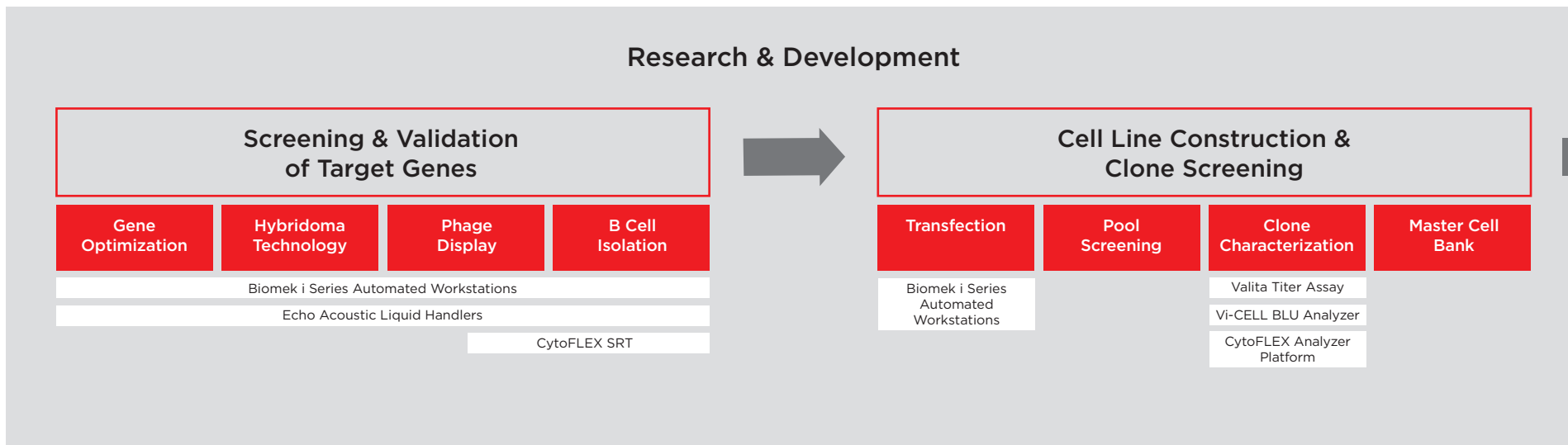
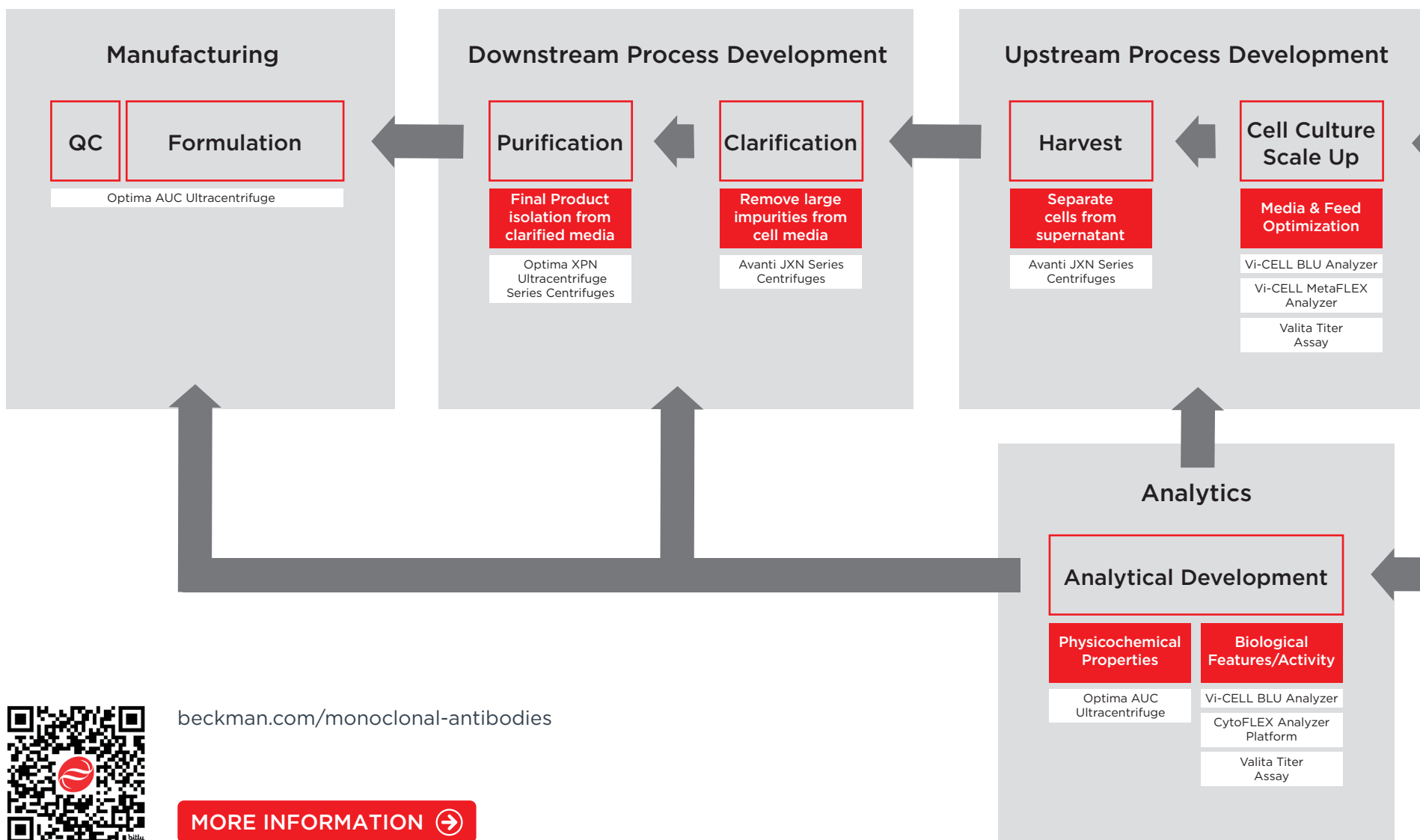


# mAbs Development Ecosystem in Biopharma

Science



Patient



**Analytics**

**Analytical Development**

Physicochemical Properties

Biological Features/Activity

Optima AUC Ultracentrifuge  
Vi-CELL BLU Analyzer  
CytoFLEX Analyzer Platform  
Valita Titer Assay



[beckman.com/monoclonal-antibodies](https://beckman.com/monoclonal-antibodies)

**MORE INFORMATION** ➔