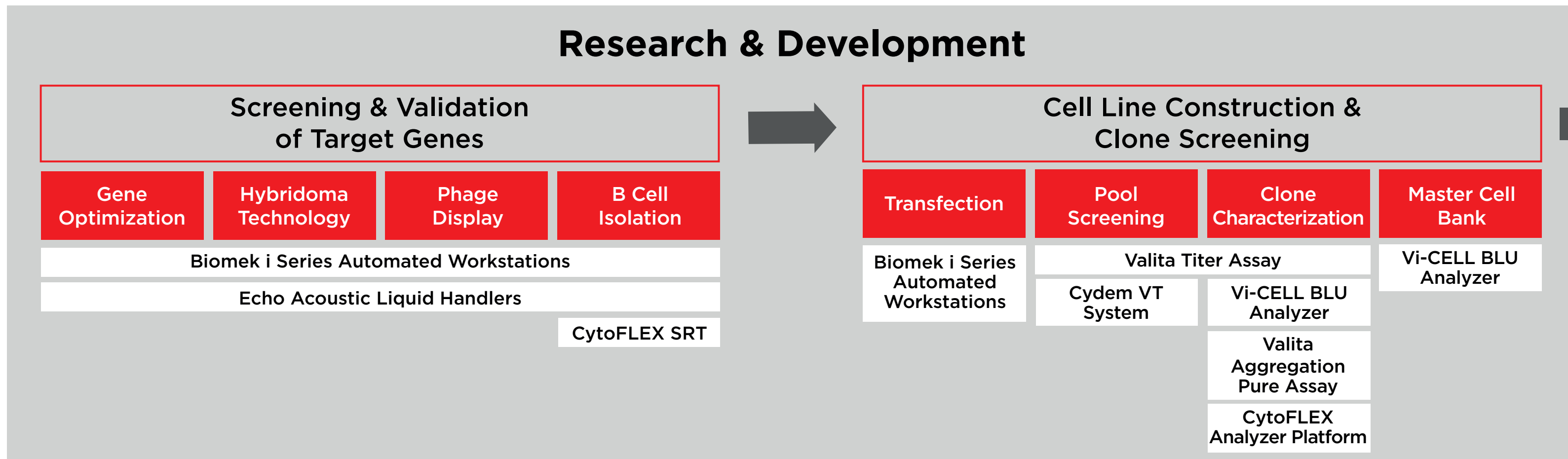


# mAbs Development Ecosystem in Biopharma

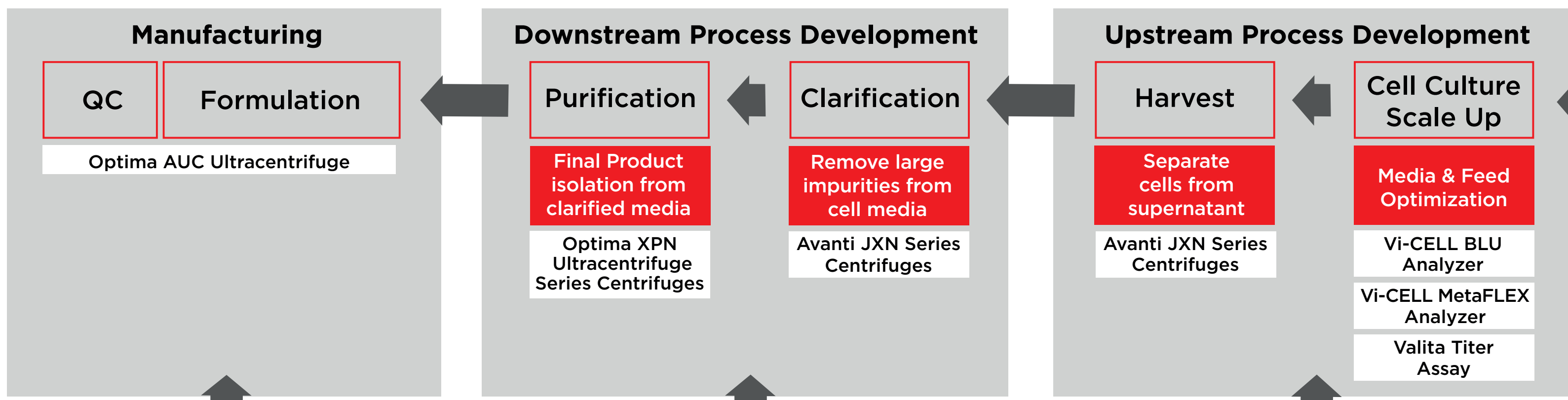
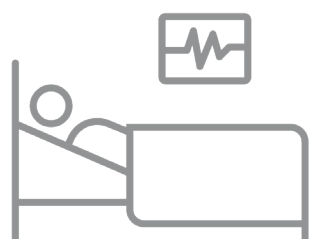
Science



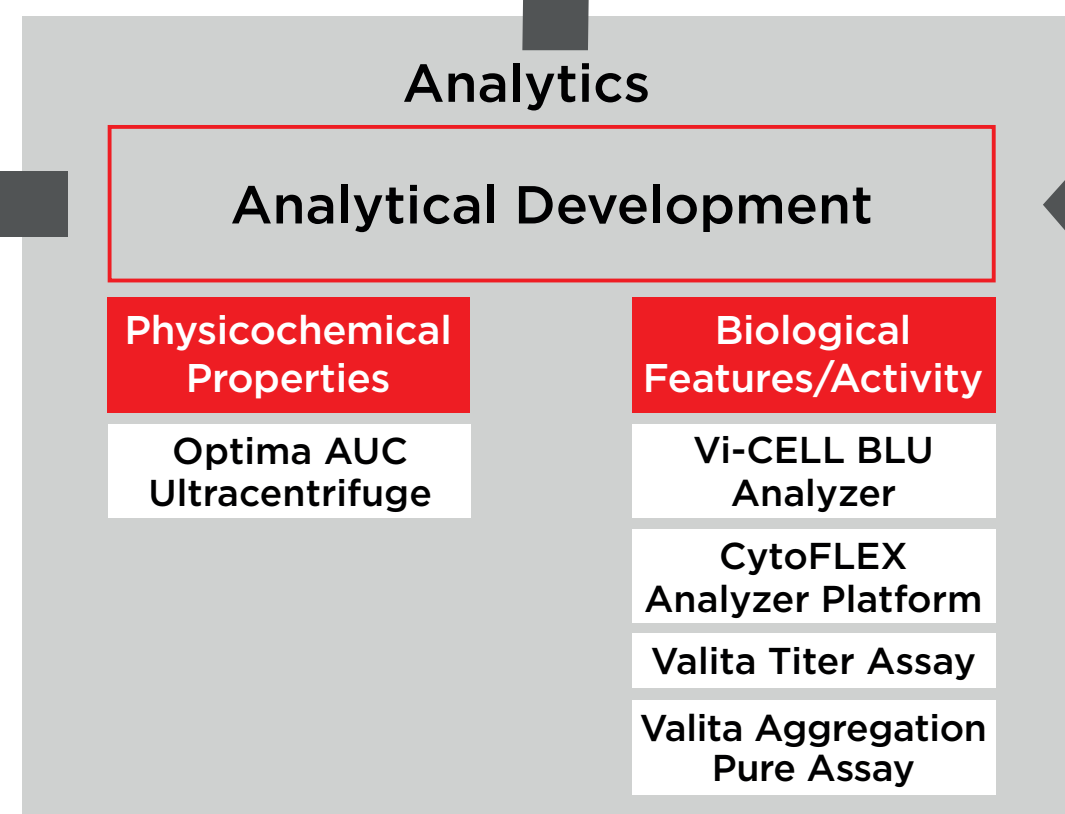
## Research & Development



Patient



## Analytics



Learn more about newest trends in the discovery, development and manufacturing of mAbs. [beckman.com/monoclonal-antibodies](https://beckman.com/monoclonal-antibodies)