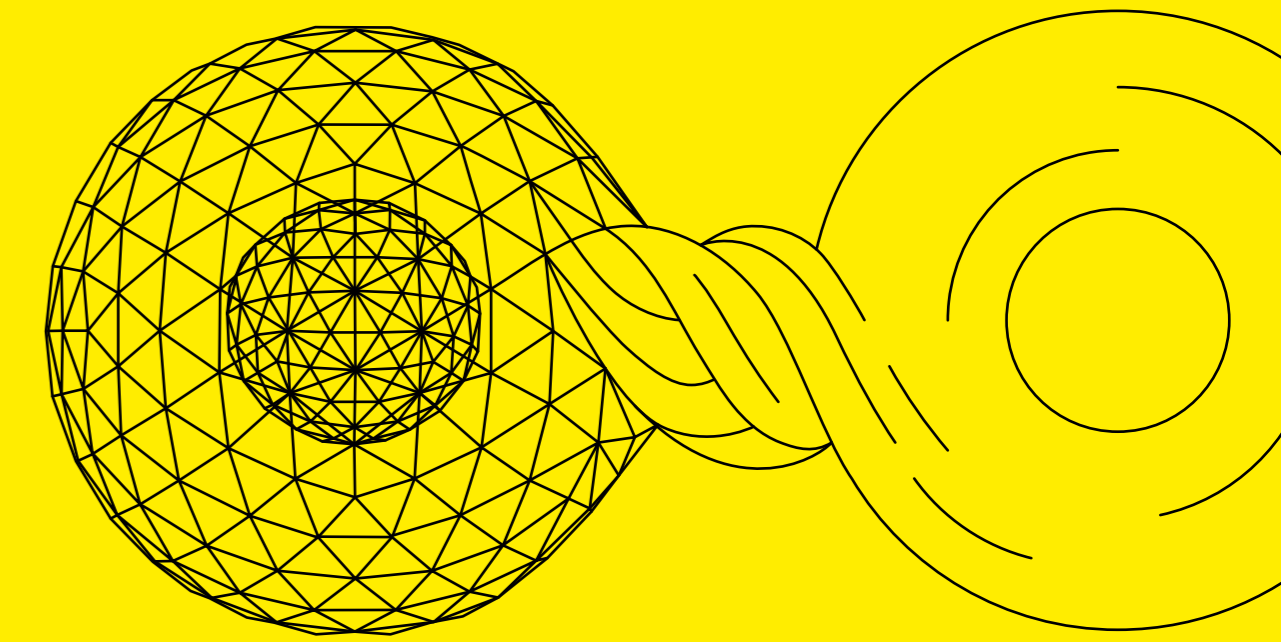


Media Formulation Optimization for Fed-Batch Processes

Decide on the Right Media up to 25% Faster
With the Help of a Virtual Bioreactor

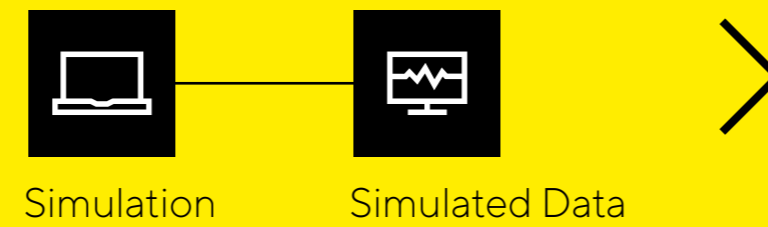


Virtual Bioreactor Media Optimization: 90 – 180 Days

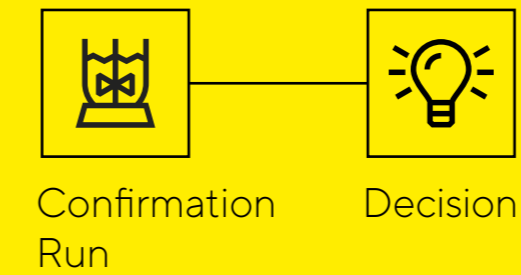
Initial Setup: 1 – 2 Months



Prediction and Analysis: 1 – 2 Months



Final: 1 – 2 Months

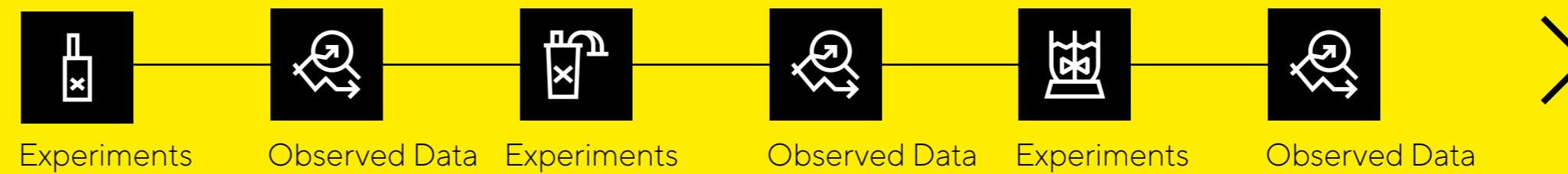


Traditional Media Optimization: 120 – 210 Days

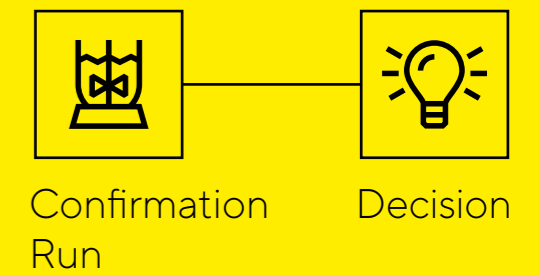
Initial Setup: 1 – 2 Months



Development Runs: 2 – 3 Months



Final: 1 – 2 Months



Illustrative Project Timeline: Theoretical project set up with the following assumptions. Initial fed-batch is 24 days from thawing first vial to harvest of first campaign. All following campaigns are 3 weeks assuming a rolling inoculum train. One week allocated for data analysis between each campaign.