

Checklist: How to Spot a Clone Worth Keeping

Use this checklist during screening and selection to confidently identify high-potential clones:

- ☐ Displays uniform morphology across all replicates and time points
- ☐ Lacks satellite colonies, doublets, or signs of contamination
- ☐ Exhibits rapid and consistent growth under standard conditions
- ☐ Produces stable and high-yield protein expression (when applicable)
- ☐ Confirmed monoclonality via white light and/or fluorescence imaging
- ☐ Selected using traceable, image-documented methods
- ☐ Falls within the optimal productivity range as defined by project criteria
- ☐ Viable for scale-up in downstream manufacturing or research processes