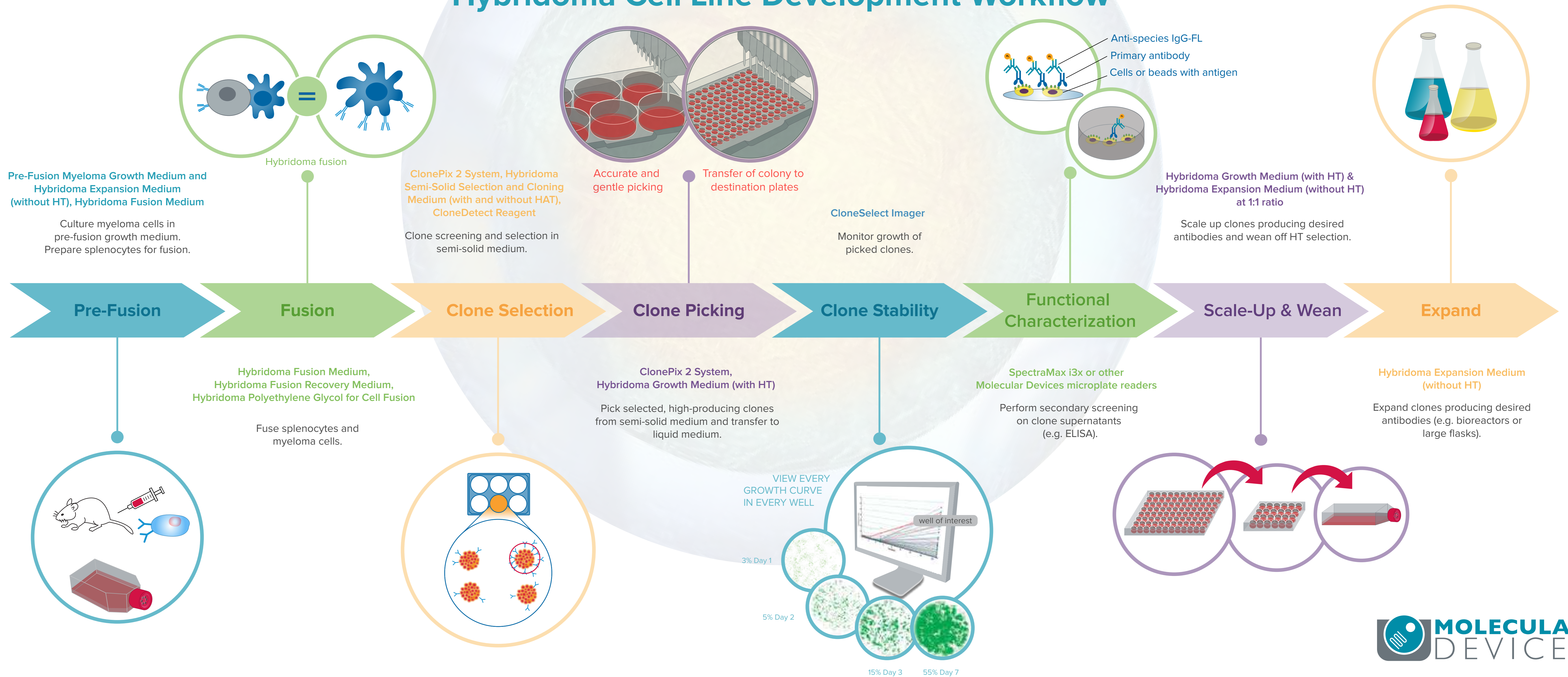


# Hybridoma Cell Line Development Workflow



To learn more, please visit [moleculardevices.com/grow-hybridomas](https://moleculardevices.com/grow-hybridomas)