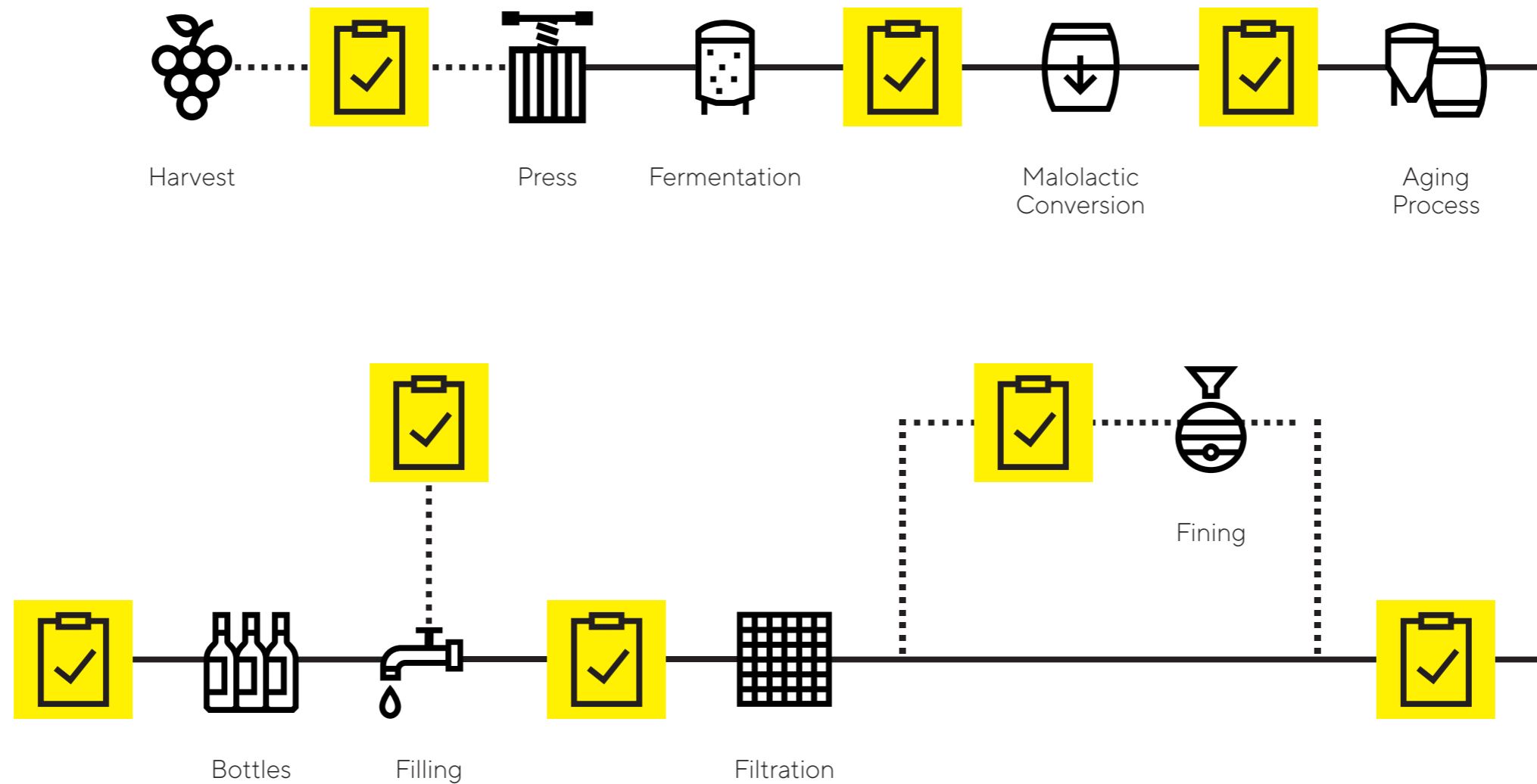


Process Flow Chart for Your Wine Process*



*This workflow may not exactly represent your winery workflow, however it generalizes the QC checkpoints.