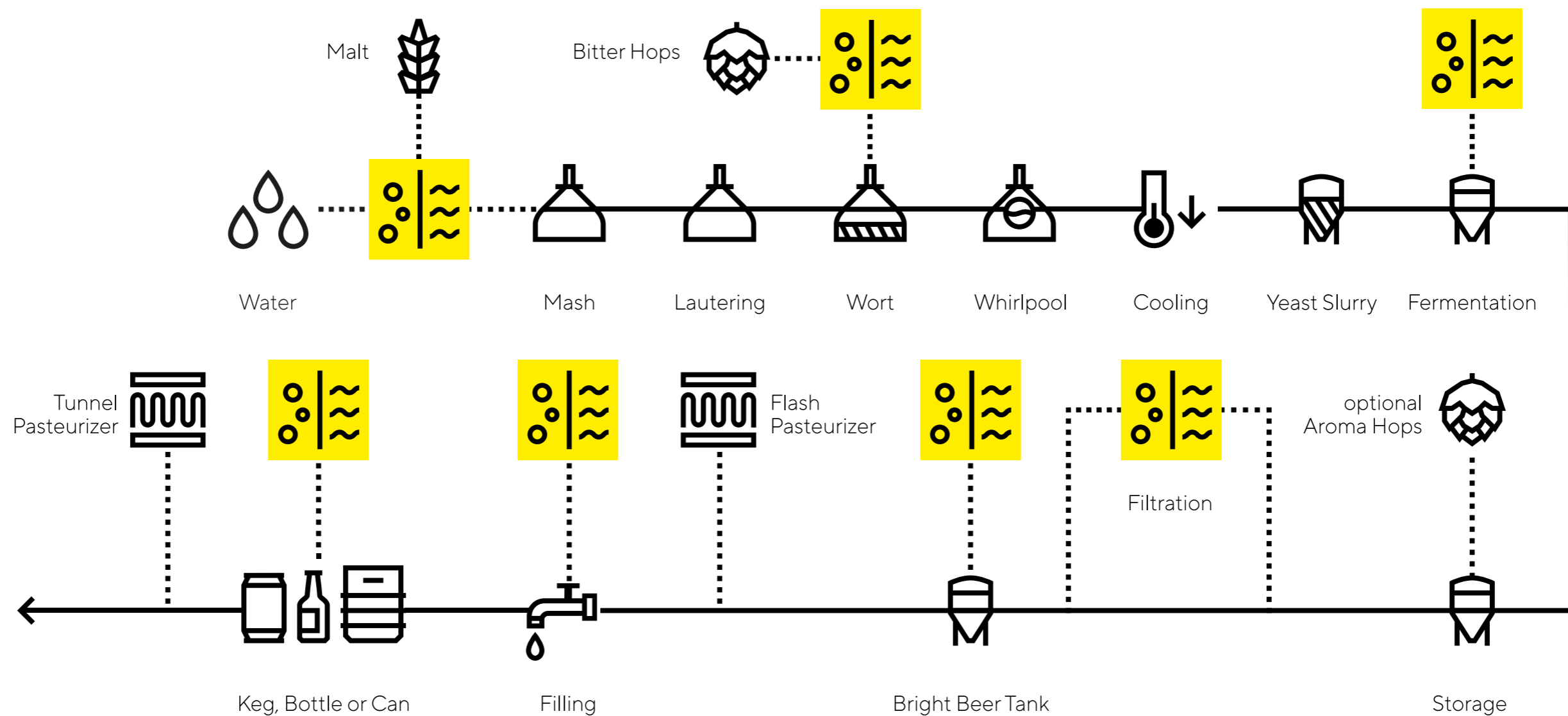


### Process Flow Chart for Your Beer Process\*



\*This workflow may not exactly represent your brewery workflow, however it generalizes the process checkpoints.