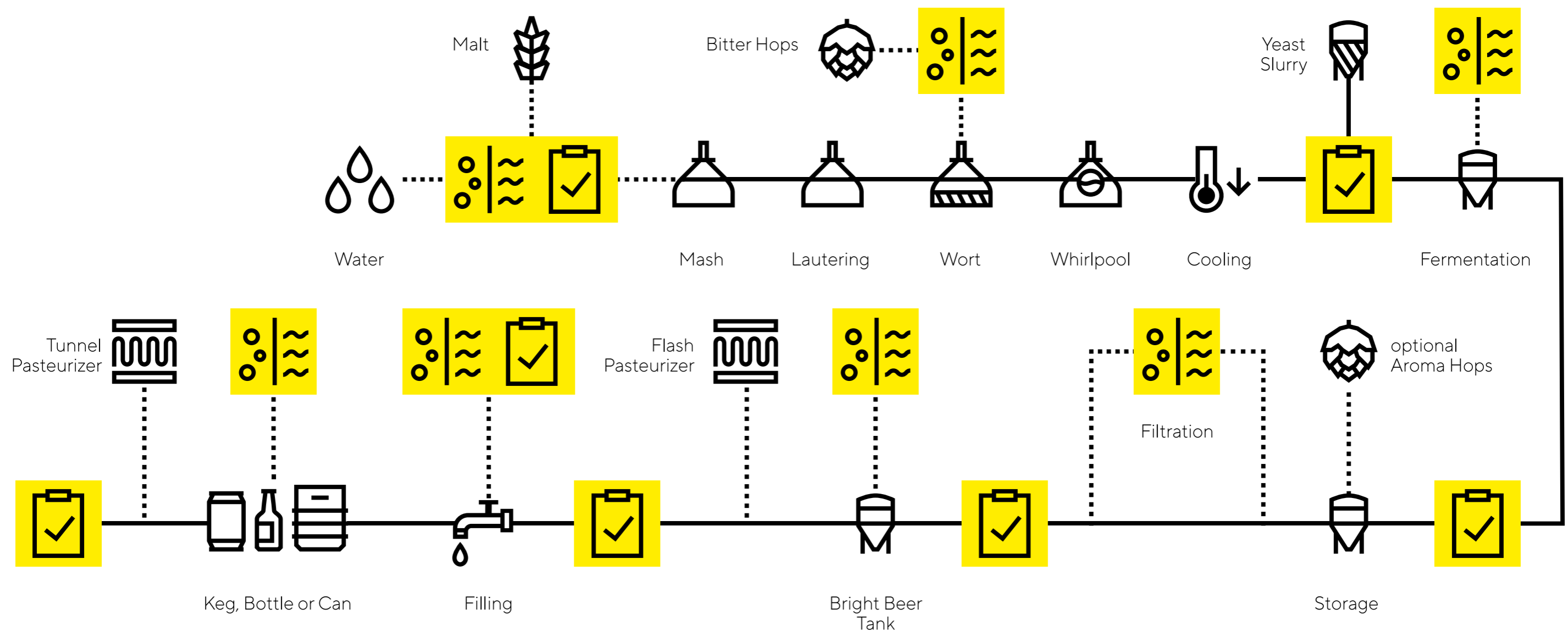


Process Flow Chart for Your Beer Process*



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