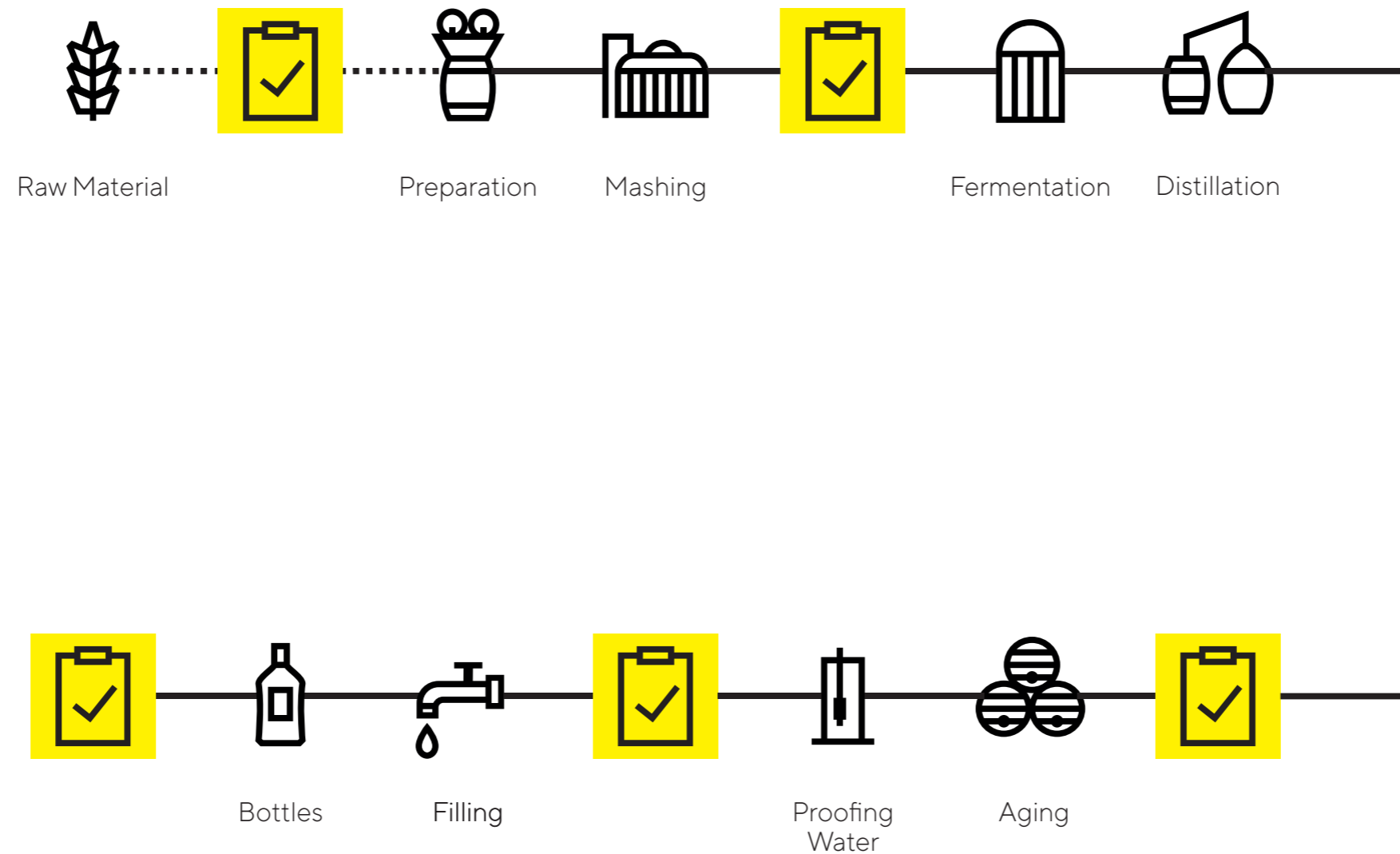


Process Flow Chart for Your Spirit Process*



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