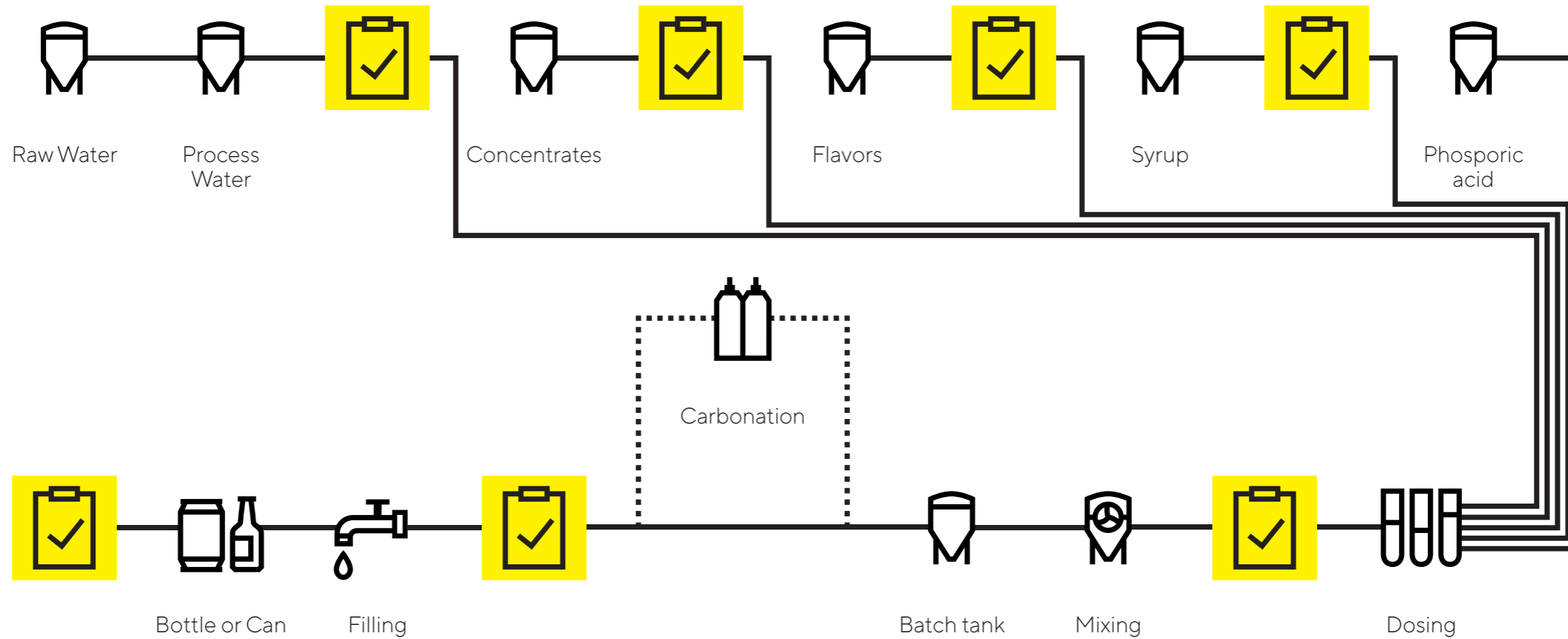


Process Flow Chart for Your Soft Drink Process*



*This workflow may not exactly represent your workflow in your soft drink manufacturing, however it generalizes the QC checkpoints.