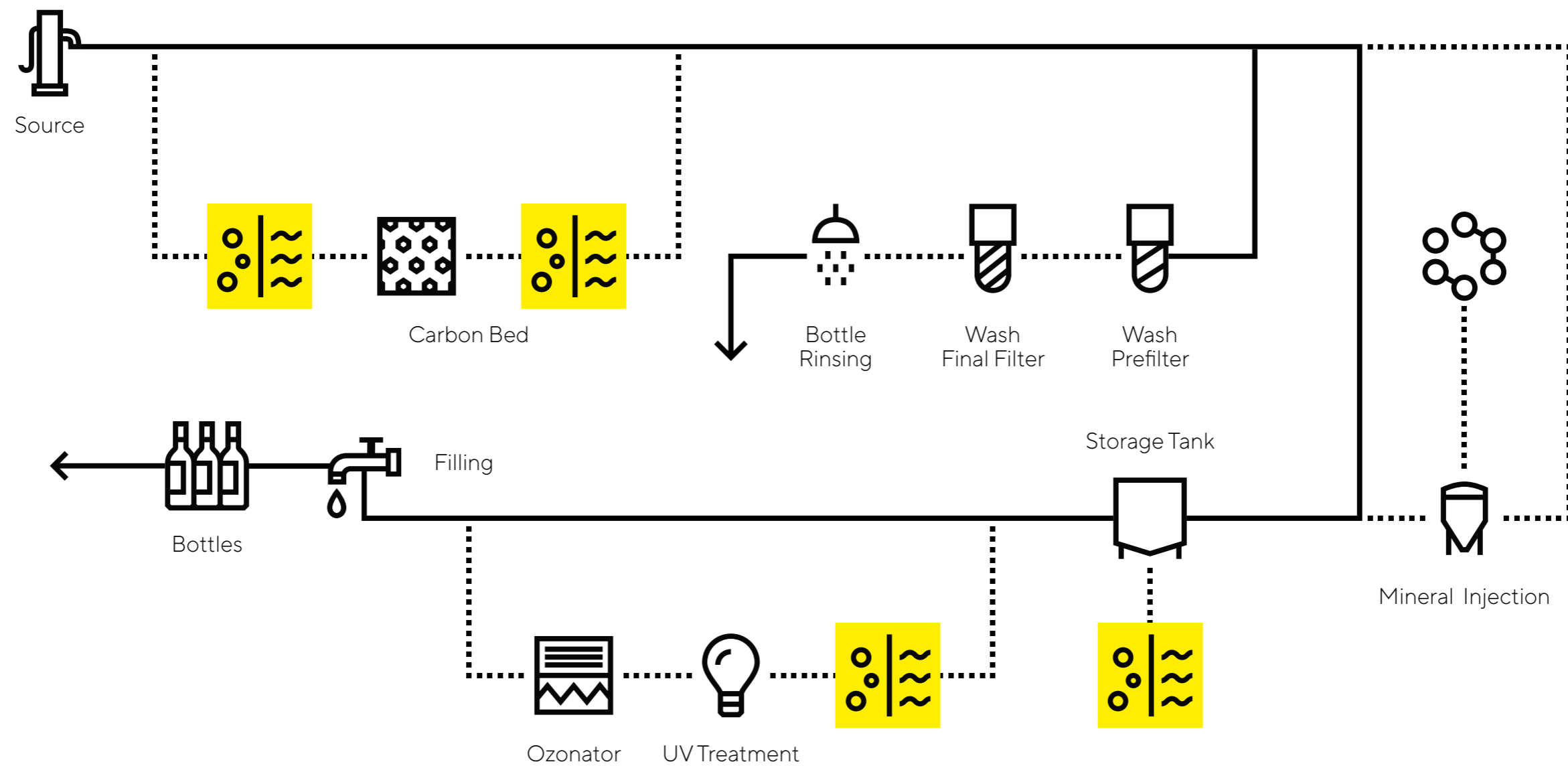


Process Flow Chart for Your Water Process*



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