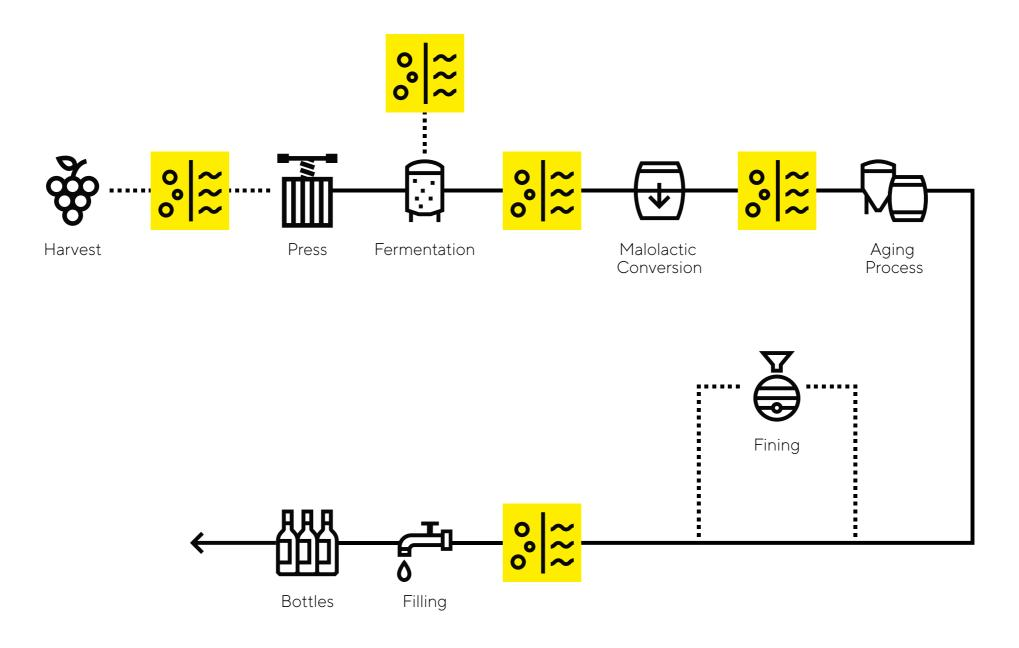
SARTORIUS

Process Flow Chart for Your Wine Process*



Simplifying Progress