

Confidence[®] Validation Service Validation Project Workflow

Your competent partner for validation

- Filter validation
- Single-use system validation
- Extractables | Leachables testing
- Virus clearance services

Quality

All validation studies follow and are compliant with regulatory requirements. We identify the right approach for each customer, resulting in over 98% initial acceptance by regulatory authorities.

Expertise

In more than 30 years Confidence[®] successfully completed over 6000 validation studies and satisfied more than 800 global customers in over 60 different countries.

Cost efficiency

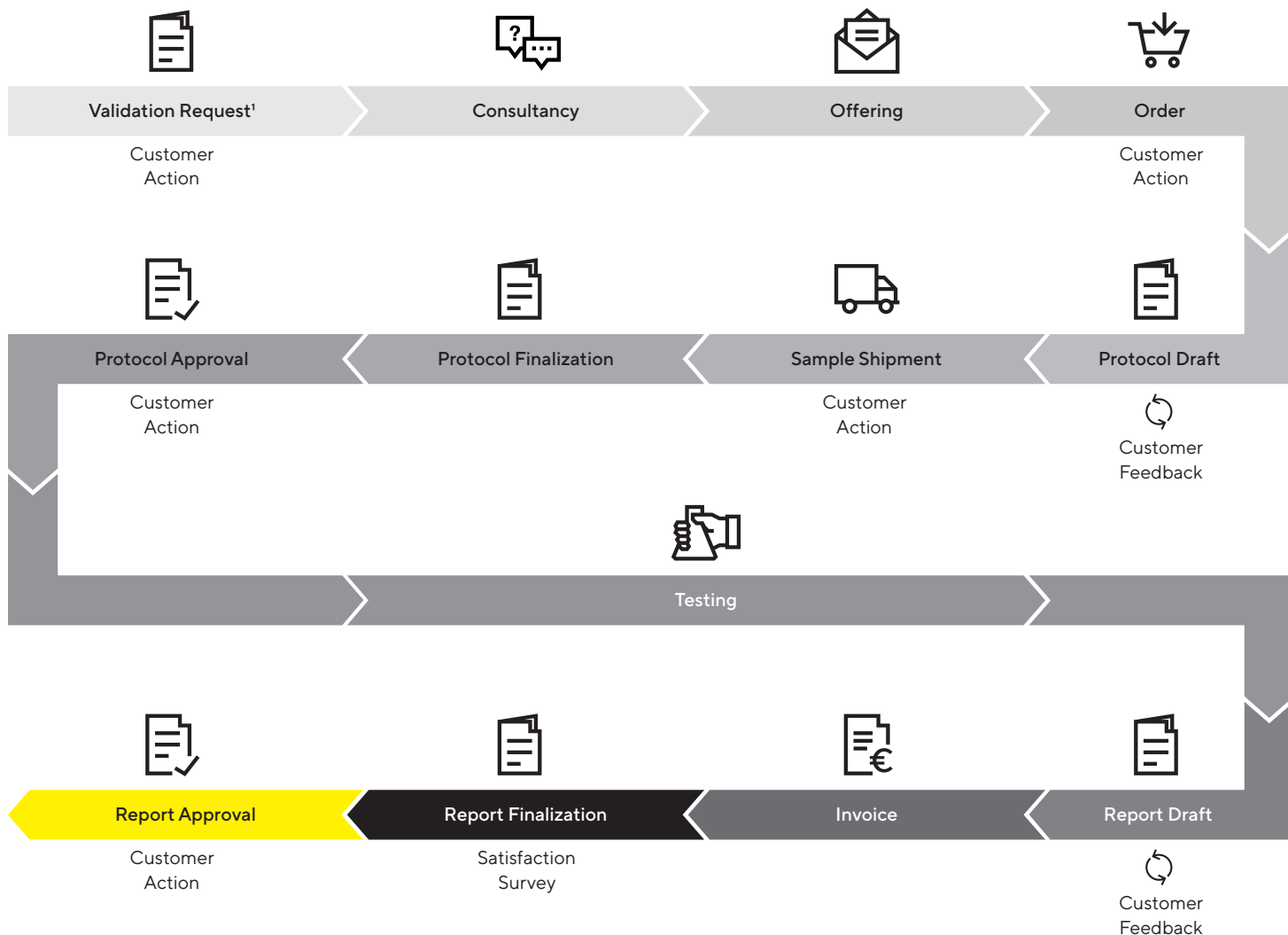
By taking a risk-based approach to each validation project, we deliver the most cost-effective testing strategy.

Fast time to market

Minimal clarification effort, agile lab scheduling and a proactive work attitude enable study times as fast as six weeks.



Confidence® Validation Services Process Chart




¹via Process Operating Parameter Sheet (POP)

Germany

Sartorius Stedim Biotech GmbH
 August-Spindler-Strasse 11
 37079 Goettingen
 Phone +49 551 308 0

USA

Sartorius Stedim North America Inc.
 565 Johnson Avenue
 Bohemia, NY 11716
 Toll-Free +1 800 368 7178

 **For more information, visit**
www.sartorius.com