



4Cell[®] CHO Platform

Cell Line Development
Service

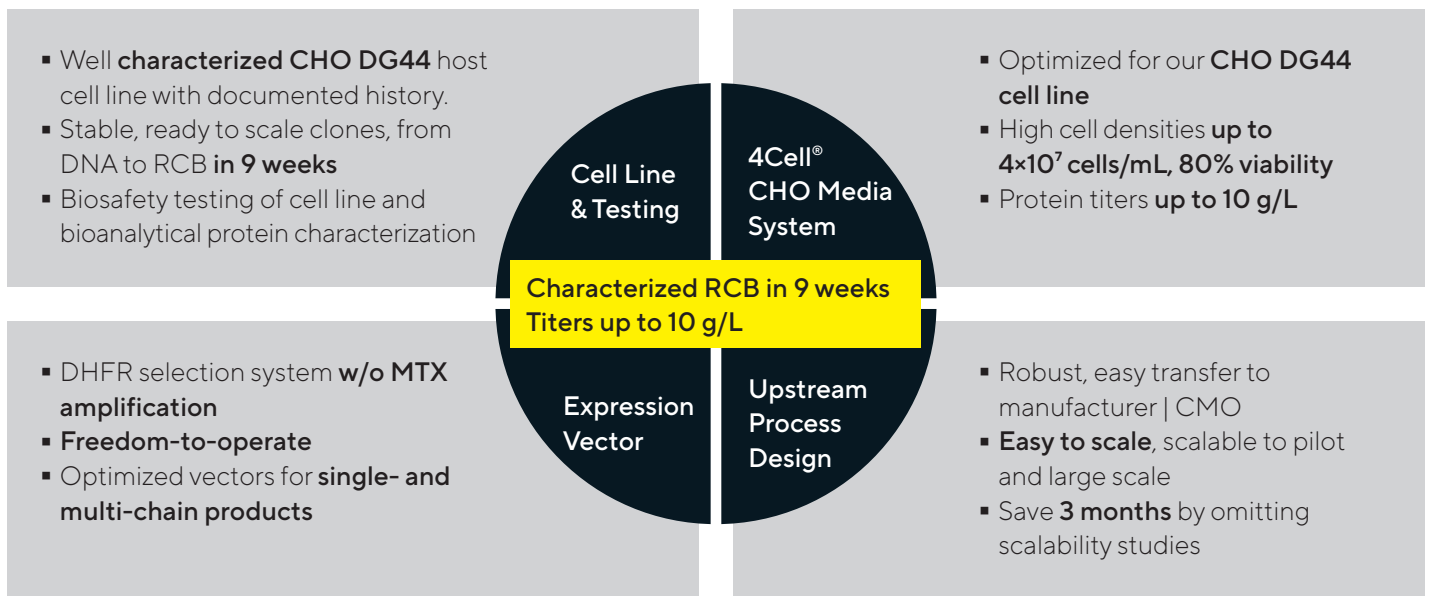
Simplifying Progress

SARTORIUS

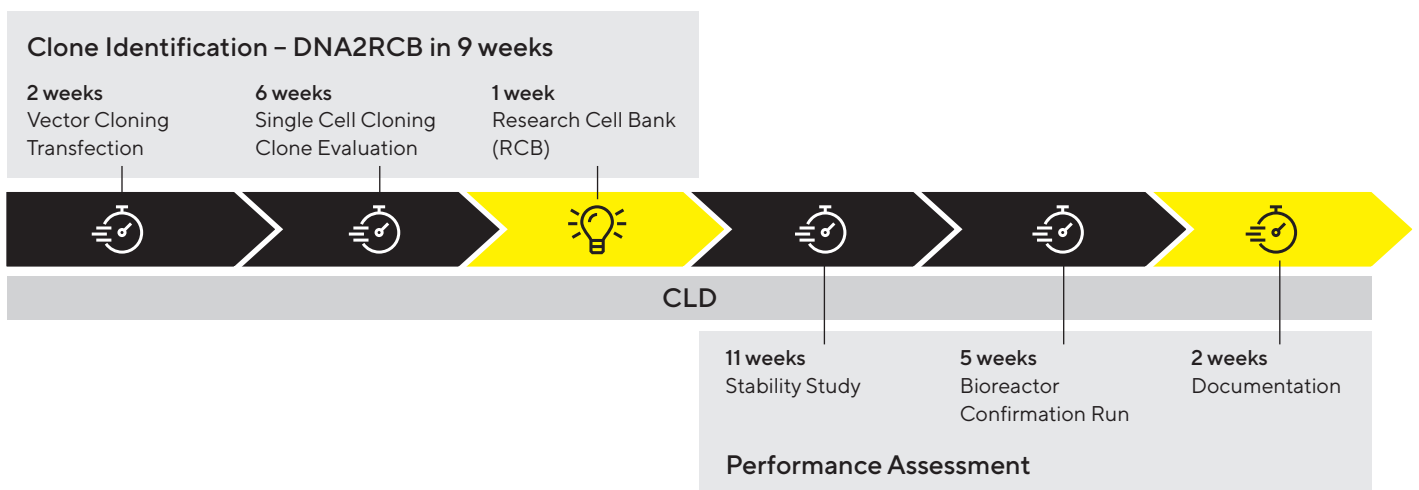
Cell Line Development Service

Sartorius is a leading provider of cell line development services allowing customers easy open access to a cost effective reliable technology platform. The Cell Line Development platform consistently delivers stable research clones from DNA to Research Cell Bank (RCB) in just 9 weeks with monoclonal antibody titers of up to 10 g/L in an easily scalable fed batch process.

Key Components



CLD: Cell Line Development, RTU: Ready to Use, RCB: Research Cell Bank, MCB: Master Cell Bank



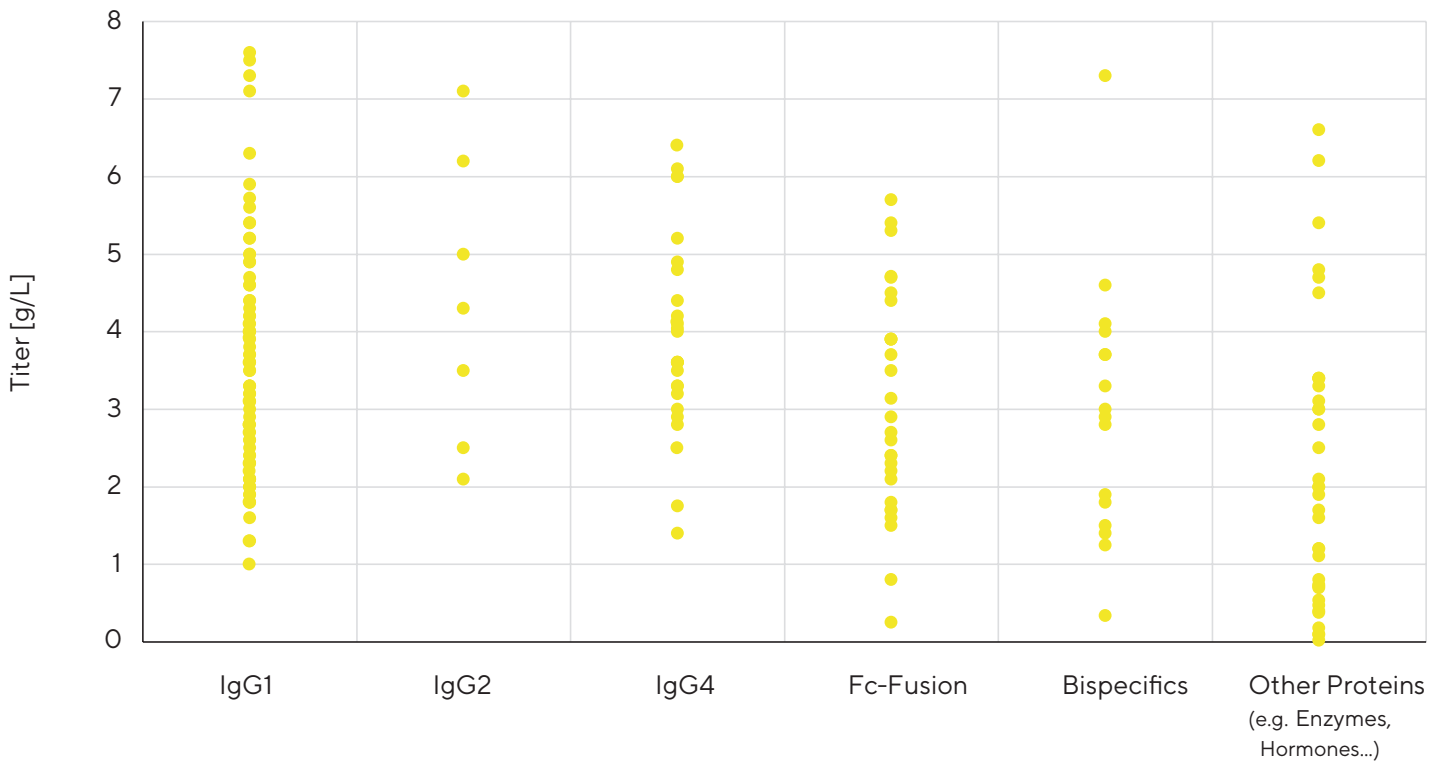
Key Benefits

Track Record and Performance

More than 240 projects using the Sartorius Cell Line Development Technology.

Track Record Cellca CLD

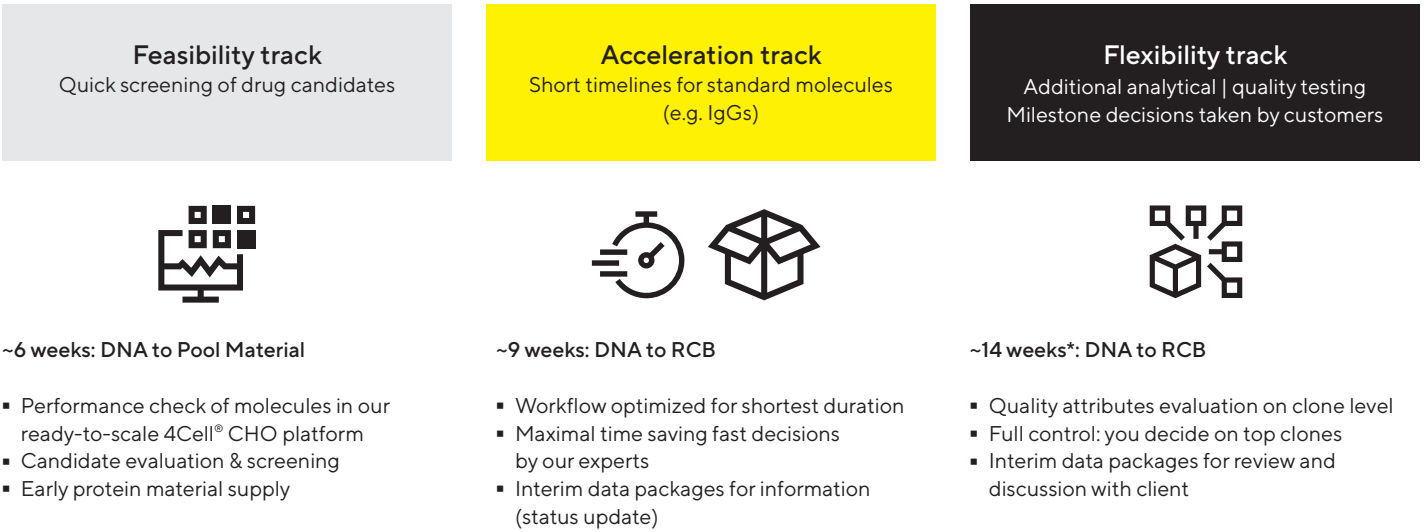
Unoptimised yields in a 12-14 fed-batch process with minimum 70% viability



Status: September 2022

Speed and Flexibility

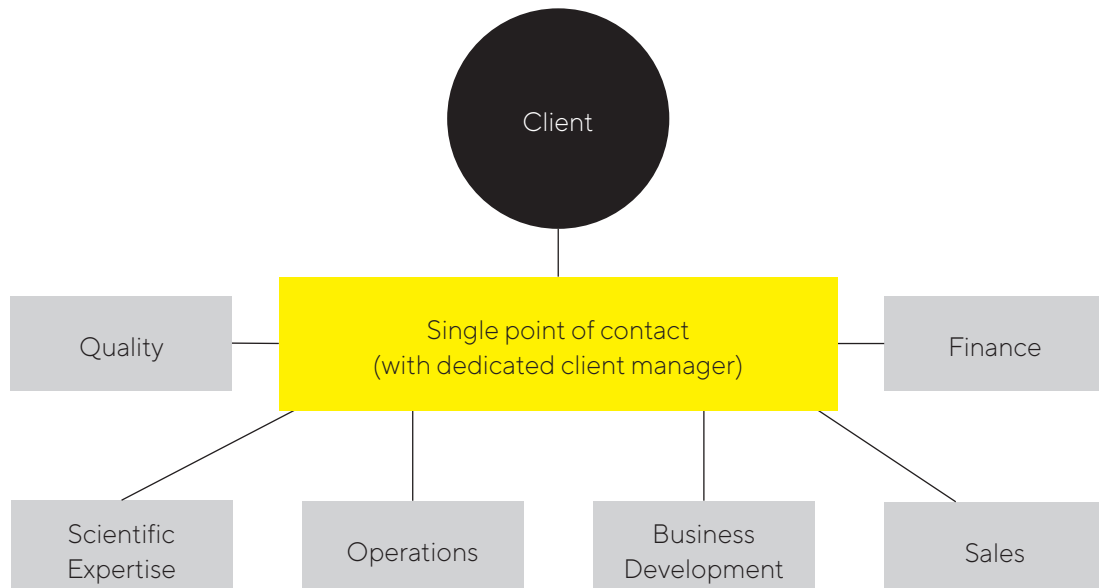
From DNA to fully characterized RCB in just 9 weeks. Save up to 3 months by omitting the need for scalability studies.



*Timeline is for a standard IgG1 NBE – scope & timeline of analytical testing might vary depending on molecule & client's needs

Customer Focus

Single point of contact throughout the whole process from DNA to fully characterized RCB.



Global Facilities of Sartorius for Cell Line Development, Cell Bank Manufacturing and Testing Services



The 4Cell® CHO Platform: Integrated Service Package from DNA to Cell Banks

DNA to Cell Banks				
Predictable performance and accelerated drug development with our integrated CHO platform				
Services	Cell Line Development	Protein Characterization	CHO Cell Banking	Biosafety testing
	<ul style="list-style-type: none"> Stable, ready to scale clones From DNA to RCB in 9 weeks Predictable protein titers: up to 10 g/L 	<ul style="list-style-type: none"> Product Quality Assessments Assay development expertise Full assay lifecycle management 	<ul style="list-style-type: none"> GMP Master & Working cell banks 500 vial @ 12 mio. cells/vial Qualified Person for cell bank release 	<ul style="list-style-type: none"> Validated GMP Release Assays Sterility, Mycoplasma, Virus Testing Genetic Characterization

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

Germany

Sartorius Stedim Cellca GmbH
Marie-Goeppert-Mayer-Str. 9
89081 Ulm-Eselsberg
Phone +49 731 206520

 For further contacts, visit
www.sartorius.com

Specifications subject to change without notice.

© 2022 Sartorius Stedim FMT S.A.S., Avenue de Jouques - Zone Industrielle Les Paluds - C.S. 91051, 13781 Aubagne Cedex, France

Publication No.: S--1539-e | Order No.: 85037-554-21 | Status: 12 | 12 | 2022