# SVISCISVS

## Product Datasheet

## Octet<sup>®</sup> Software Version 13

## Added Features to Improve EC<sub>50</sub>, IC<sub>50</sub> and Potency Analysis

### New Key Features

- New EC<sub>50</sub> and IC<sub>50</sub> analysis capabilities to determine drug product activities using surrogate potency assays, reducing assay variability, timelines, and costs
- Data bridge to support data transfer between Sartorius Octet<sup>®</sup> and Sartorius Ambr<sup>®</sup> bioreactor systems to accelerate cell line development workflows

#### Overview

Sartorius Octet<sup>®</sup> BLI Discovery and Analysis Studio Software are evolving to address growing applications, and user needs with every release version.

## Dose-Response Analysis Capabilities

Determination of product activity is vital during drug development and manufacturing to ascertain the intended biological activity of the therapeutic. A potency or potencyindicating assay is a requirement by regulatory bodies before the release of every manufactured drug lot. Typically, potency measurements are obtained using bioassays in animal models or cell-based assays to capture biological activities directly. However, alternative assay formats such as ELISA or BLI are used for potency measurements when drug binding activities to target molecules correlate with the biological activity of the drug. These assays are coined as potency-indicating or surrogate assays for their ability to reveal drug activity based on target binding. Octet® doseresponse assays provide fast time to results with automated workflows, minimal hands-on times, and reduced assay variabilities. Unlike ELISA, all assay steps can be monitored and evaluated in real time.

- Conveniently setup EC<sub>50</sub> and IC<sub>50</sub> experiments with built-in experimental method templates in Octet<sup>®</sup> Discovery Software.
- 3-PL, 4-PL, 5-PL, linear, and semi-log fitting algorithms
- Monitor biosensor loading responses and flag data points based on user-defined biosensor loading criteria
- Independent and global dose-response fitting options and similarity assessments to effectively compare against standard and test sample groups
- Create customized data reports for presenting and reporting dose-response data and conclusions

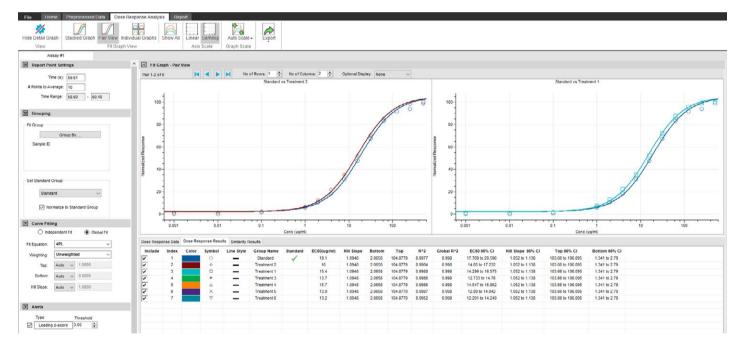


Figure 1: Anti-Her2 lot-to-lot potency evaluation using an Octet<sup>®</sup> potency assay. Relative potency evaluation between two anti-Her2 antibody lots using the binding activity to Her2. Dose-repose binding data from multiple lots were fitted globally to a 4-PL fitting algorithm in Octet<sup>®</sup> Analysis Studio Software.

## Octet<sup>®</sup>- Ambr<sup>®</sup> 15 Data Bridge

Octet<sup>®</sup> BLI platforms are routinely implemented in upstream and downstream cell line development workflows measuring critical quality attributes (CQAs) such as product titer, binding activity, or glycosylation ranking. The Ambr<sup>®</sup> 15 is an automated high-throughput single-use bioreactor system that provides parallel operation of up to 48 single-use microbioreactors at a time where multiple cell culture experiments can be used to evaluate different cell lines, clones, or to identify desirable cell culture process parameters. Ambr<sup>®</sup> cell culture workflows are often combined with label-free binding assays due to fast time to results and minimal hands-on time.

- Built-in Octet<sup>®</sup>-Ambr<sup>®</sup> data bridge in Octet<sup>®</sup> BLI Discovery Software for importing Ambr<sup>®</sup> sample and condition information, saving analyst time and eliminating data transfer errors.\*
- Directly import Ambr® 15 sample plate map configurations with experimental information, rearrange sample well positions for Octet® assays, and modify sample information or replicate groups using the Octet®-Ambr® 15 data bridge.
- Preconfigure data import and export file locations to streamline data import and export to and from the Octet<sup>®</sup> to Ambr<sup>®</sup> 15 data bridge.
- Seamless transfer of Octet<sup>®</sup> data analytics from Octet<sup>®</sup> Analysis Studio Software to Ambr<sup>®</sup> Software for integrating cell culture analytics for faster data feedback.\*

Ambr® Data Import							1 Rate Definition 2 Sensor Assignment 3 Review Experiment 4 Run Experiment			
sex Impo	t Original Well	ell Plate	Station V	essel Timestamp System ID	Experiment B	A V 1 2 3 4 5 6 7 8 9 10 11 12	in this step, all the information about the sample plate and its v First, check the assay settings. Then highlight one or more we	els will be entered. Is on the sample plate, and right-click to enter.	modfy well data.	
	A1 .	Sample	ate CS1+2 CS1 1	05/May/2022 07:34 ambr48-R5-4	20220506_TDT61-PD-AR8	ADOOOOOOOOOOOOOOOO	Read Head Schannels high sensitivity)	Plate 1 Table (96 wells)		
	A2	Sample	ate CS1+2 CS1 2	09/May/2022 07:36 ambr48-R5-4	2 20220506_TDT61-PD-AR8	BOODOOOOOO		Concentration units: up liel 🗸	Export. Import. Print.	
	A3	Sample	ate CS1+2 CS1 3	05/May/2022 07:37 ambr48-R5-4	2 20220506_TDT61-PD-AR8		Acquistion Rate Standard (5.0 Hz) V	Well Sample ID Replicate Group	Type Conc (pg/ml) Dilutio	on Factor Information
	A4		ate CS1+2 CS1 4	09/May/2022 07:38 ambr48-R5-4			Assay Settings Assay: Basic Quantitation with Regeneration Modely	C3 C51-1	Ukknown n/a	anbr48-R5-452.20220506_TDT61-PD-AR8.CS1.1.10/May/202
	AS		ate CS1+2 CS1 5	09/May/2022 07:39 ambr48-R5-4			Single analy Single analy Countration Representation Representation Single analy Time (b) Representation Representation Single analy Representation Single analy Representation Single analy Representation Representation Single analy Representation	D3 CS1-7	Utiknown n/a	anbr48/R5-452.20220506_TDT61-PD-AR8,CS1.7.10/May/202
								E3 CS1-1	O Unknown n/a	anbr48-R5-452.20220506_TDT61-PD-AR8.C51.1.11/May/202
· 🖓	AG		ate CS1+2 CS1 6	09/May/2022 07:40 ambr48-R5-4				F3 CS1-7	Utiknown n/a	anbr45/R5-452.20220506_TDT61/PD-AR8.CS1.7.11/May/202
	81	Sample	ate CS1+2 CS1 7	09/May/2022 07:41 ambr48-R5-4	20220506_TDT61-PD-AR8	FOODOOOOO		A4 CS1-2	Usknown n/a	ambr48/R5-452.20220506_TDT61/PD AR8.CS1.2.09/May/202
	82	Sample	ate CS1+2 CS1 8	09/May/2022 07:42 ambr48-R5-4	2 20220506_TDT61-PD-AR8			84 CS1-8	Utknown n/a	anbr45/R5-452,20220506_TDT61-PD-AR8.C51.8.09/May/20
R	83		ate CS1+2 CS1 9	09/May/2022 07:43 ambr48-R5-4		GOOOOOOOOOOOO	3 cycles per regeneration Pre-conditioning Enabled	C4 C51-2	Utknown n/a	ambr48/R5-452,20220506_TDT61/PD-AR8.CS1,2:10/May/20
	84		ate CS1+2 CS1 10	09/May/2022 07:44 ambr48-R5-4		100000000000000		D4 C51-8	<ul> <li>Usknown n/a</li> </ul>	ambr48-R5-452,20220506_TDT61-PD-AR8.C51.8.10/May/20
						<b>1</b> 000000000000000000000000000000000000	Plate 1 (95 wells) V Modily	E4 CS1-2	Ukknown n/a	anbr45-R5-452.20220506_T0T61-PD-AR8.C51.2.11/May/2
	85		ate CS1+2 CS1 11	09/May/2022 07.45 ambr48-R5-4			V 1 2 3 4 5 6 7 8 9 10 11 12	F4 CS1-8	<ul> <li>Unknown n/a</li> </ul>	ambr48/R5-452,20220506_TDT61/PD-AR8,CS1.8,11/May/2
	86	Sample	ate CS1+2 CS1 12	2 09/May/2022 07:47 ambr48-R5-4	20220506_TDT61-PD-AR8			A5 CS1-3	Utknown n/a	anbr48-R5-452,20220506_TDT61-PD-AR8,CS1,3,09/May/20
	C1 (	Sample	ate CS1+2 CS1 1	10/May/2022 07:34 ambr48-R5-4	2 20220506 TDT61-PD-AR8			85 CS1-9	Usknown n/a	ambr48-R5-452,20220506_TDT61-PD-AR8.C51.9.09/May/20
	62	Sample	ate CS1+2 CS1 2	10/May/2022 07:36 ambr48-R5-4	2 20220506_TDT61-PD-AR8	Ready to import		C5 C51-3	<ul> <li>Unknown n/a</li> </ul>	anbr48/R5-452,20220506_TDT61/PD-AR8.CS1.3.10/May/20
	3		ate CS1+2 CS1 2 ate CS1+2 CS1 3	10/May/2022 07:37 ambr48-R5-4				05 CS1-9	Ukknown n/a	anbr48-R5-452.20220506_TDT61-PD-AR8.C51.9.10/May/20
1.4								E5 CS1-3	Utknown n/a	ambr48/R5-452,20220506_TDT61/PD-AR8.CS1.3.11/May/20
	C4 (	Sample	ate CS1+2 CS1 4	10/May/2022 07:38 ambr48-R5-4	2 20220506_TDT61-PD-AR8			F5 C51-9	Unknown n/a	ambr48-R5-452,20220506_TDT61-PD-AR8.C51.9.11/May/20
	C5 (	Sample	ate CS1+2 CS1 5	10/May/2022 07:39 ambr48-R5-4	2 20220506_TDT61-PD-AR8			A6 CS1-4	Unknown n./a	ambr48-R5-452,20220506_TDT61-PD-AR8.C51,4.09/May/20
R	C6 (	Sample	ate CS1+2 CS1 6	10/May/2022 07:40 ambr48-R5-4	2 20220506_TDT61-PD-AR8			B6 CS1-10	Utiknown n/a	anbr45/R5-452,20220506_TDT61/PD-AR8.C51.10.09/May/2
	D1		ate CS1+2 CS1 7	10/May/2022 07:41 ambr48-R5-4				C6 C514	Utknown n/a	anbr45-R5-452,20220506_TDT61-PD-AR8,CS1,4.10/May/20
				10/May/2022 07:42 ambr48-R5-4				D6 CS1-10	Utknown n/a	anbr48-R5-452.20220506_TDT61-PD-AR8.CS1.10.10/May/2
10 🔽 D2	02	Sample	Sample plate CS1+2 CS1 8	rurmay/2022/07/42/ambr40-H5-4	20220506_101614P0-AH8	£		E6 C51-4	Unknown n/a	anbr45/R5-452,20220506_TDT61/PD-AR8,C51,4,11/May/20
					>		Standard O Control Unassigned	F6 CS1-10	O Unknown n/a	anbr45-R5-452.20220506_TDT61-PD-AR8.CS1.10.11/May/2
							O Unknown O Reserved	A7 C51-5	O Urknown n/a	anbr45-R5-452.20220506_TDT61-PD-AR8.C51.5.09/May/20
Choose 8	les				Clear			87 CS1-11 C7 CS1-5	O Unknown n/a	ambr48 R5452 20220506_TDT61 PD AR8.CS1.11.09/May/2
			100					C7 C51-5 D7 C51-11	O Unknown n/a	ambr48-R5-452.20220506_TDT61-PD-AR8.C51.5.10/May/20
Information mapping			Well mapping					O Utknown n/a	anbr48/R5-452.20220506_TDT61/PD-AR8.C51.11.10/May/2	
				on the import table to apply a well mapping	Ue			E7 CS1-5	O Urknown n/a	anbr48/R5-452.20220506_TDT61/PD AR8.CS1.5.11/May/20
ple	Culture Static	Vessel (CS1-1)	24 1	vell to 96				F7 C51-11	O Urknown n/a	ambr48-R5-452.20220506_TDT61-PD-AR8.CS1.11.11/May/
Acate	deave blank)		Direct	IL TR	Left Fight			A8 C516	Utknown n/a	anbr48/R5-452.20220506_TDT61/PD-AR8.CS1.6.09/May/2
	tergine (reality)				tion in the local division of the local divi			B8 C51-12	Utknown n/a	anbr45-R5-452.20220506_TDT61-PD-AR8.C51.12.09/May/2
				BL BR	Down	Import		C8 C51-6 08 C51-12	O Unknown n/a	ambr45/R5-452.20220506_T0T61/PD-AR8.C51.6.10/May/ ambr45/R5-452.20220506_T0T61/PD-AR8.C51.12.10/May/

Figure 2: Ambr® 15 process data import to Octet® BLI Discovery Software for product concentration measurement in the Octet® system. (A) The new data interface allows for importing critical sample information from Ambr® 15 process runs. Users can track sample information and reconfigure sample plate maps suitable for the Octet® assay without manually entering information in the Octet® Software. (B) Ambr® 15 sample information is automatically populated for an Octet® experiment and method set up in Octet® BLI Discovery Software.

#### New Virus Concentration Units

Setup, calculate, and report virus concentration data in VP/mL, VG/mL, CP/mL and GC/mL units in Octet® BLI Discovery and Analysis Studio Software.

### Reduce Experiment Downtimes Even Further on Octet® RH96 Instruments

Octet<sup>®</sup> RH96 instruments are powered by multiple lamps supporting the 96-channel read heads. In the event of a malfunctioning lamp, the Octet<sup>®</sup> BLI Discovery Software will guide users to continue using the instrument with the remaining lamps and reduced throughput until service on the Octet<sup>®</sup> instrument is completed to reduce instrument downtimes.

### Specifications

Component	Requirement
Processor	3.0 GHz Quad Core processor or greater
Operating system	Windows 10 or Windows 11 Professional, 64-bit
Memory	4 GB RAM or greater
Graphic display	Monitor with 1920x1080 resolution or better
Hard disk	250 GB hard drive or greater. 200 MB available hard disk space required for application files more to acquire data – a single experiment can range from 1 MB to over 100 MB.
Additional software	Microsoft Office

\*Use of Octet<sup>®</sup>-Ambr<sup>®</sup> 15 Data Bridge will require purchase of Ambr<sup>®</sup> 15 data interface license and require Ambr<sup>®</sup> Software version R10.0 or above.

## Ordering Information

Part No.	Description			
50-5029	Upgrade of Octet® instrument software to the latest version of software. Includes 4 seat licenses. Includes Octet® Software Version 13. Requires previous version of Octet® Software.			
50-5098	4-seat licenses of Octet® Software version 13.X. Includes FREE upgrades within 13.X. Requires previous purchase (50-5029) or installation of version 13.X software.			
50-5099	1-seat license of Octet® Software version 13.X. Includes FREE upgrades within 13.X. Requires previous purchase (50-50 or installation of version 13.X software.			
50-5100	2-seat license of Octet® Software version 13.X. Includes FREE upgrades within 13.X. Requires previous purchase (50-5029) or installation of version 13.X software.			
50-5101	Site license for Octet® Software version 13.X for offline data analysis. Installation and use of the Octet® Software vers 13.X at a physical site (single city or ZIP code) as defined in the purchase order. Includes FREE upgrades within 13.X. Requires previous purchase (50-5029) or installation of version 13.X software.			
50-5034	Upgrade of Octet <sup>®</sup> CFR instrument software to the latest version of software. Includes 4 seat licenses. Requires previou version of Octet <sup>®</sup> CFR Part 11 Software.			
50-5102	4-seat licenses of Octet® BLI CFR part 11 software version 13.x at a physical site (single city or ZIP code) as defined in purchase order. Includes FREE upgrades within 13.X. Requires previous purchase of Octet® BLI CFR part 11 Software version 13.x.			
50-5103	1-seat licenses of Octet® BLI CFR part 11 Software version 13.x at a physical site (single city or ZIP code) as defined in the purchase order. Includes FREE upgrades within 13.X. Requires previous purchase of Octet® BLI CFR part 11 Software version 13.x.			
50-5104	2-seat licenses of Octet® BLI CFR part 11 Software version 13.x at a physical site (single city or ZIP code) as defined i purchase order. Includes FREE upgrades within 13.X. Requires previous purchase of Octet® BLI CFR part 11 Softwa version 13.x or previous installation of a non-CFR Octet® software. Includes 2 licenses.			
50-5105	Site license for Octet® BLI CFR part 11 Software version 13.x at a physical site (single city or ZIP code) as defined in the purchase order. Includes FREE upgrades within 13.X. Requires previous purchase of Octet® BLI CFR part 11 Software version 13.x.			
50-0300	13.X SOFTWARE VALIDATION PACKAGE. For researchers working in GLP or GMP laboratories, the Octet® Software Validation Package provides comprehensive documentation and tools available to validate GxP administrator features, and quantitation and kinetics functions for data generated on Octet® instrumentation.			
001-8B39	Ambr® 15 Octet® data interface.			

#### Germany

USA

Sartorius Lab Instruments GmbH & Co. KG Otto-Brenner-Strasse 20 37079 Goettingen Phone +49 551 308 0 Sartorius Corporation 565 Johnson Avenue Bohemia, NY 11716 Phone +1 888 OCTET 75 Or +1 650 322 1360

For further contacts, visit www.sartorius.com/octet-support