

**Biomek i7 Dual
Hybrid Workstation
Automated Cytiva™
Sera-Xtracta
Virus/Pathogen Kit**



2023-GBL-EN-100795-v1

Disclaimer

Biomek Automated Workstations are not intended or validated for use in the diagnosis of disease or other conditions.

The Cytiva™ Sera-Xtracta Virus/Pathogen kit is intended for research use only. It is not intended for use in any clinical or *in vitro* procedures for diagnostic purposes. Do not use internally or externally in humans or animals.

© 2023 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries. All other trademarks are the property of their respective owners.

Beckman Coulter makes no warranties of any kind whatsoever express or implied, with respect to this protocol, including but not limited to warranties of fitness for a particular purpose or merchantability or that the protocol is non-infringing. All warranties are expressly disclaimed. Your use of the method is solely at your own risk, without recourse to Beckman Coulter. This protocol is for demonstration only; and is not validated by Beckman Coulter.

Cytiva™ Sera-Xtracta Virus/Pathogen Kit

- The Cytiva™ Sera-Xtracta Virus/Pathogen kit uses bead-based technology to isolate viral/bacterial nucleic acid (RNA and DNA).
- It is optimized to isolate nucleic acid from swab samples collected in universal transport media.
- 100µL-400µL of transport media can be purified per sample.
- Superparamagnetic SeraSil-Mag™ beads selectively bind total nucleic acid while removing unwanted material efficiently through quick wash steps.

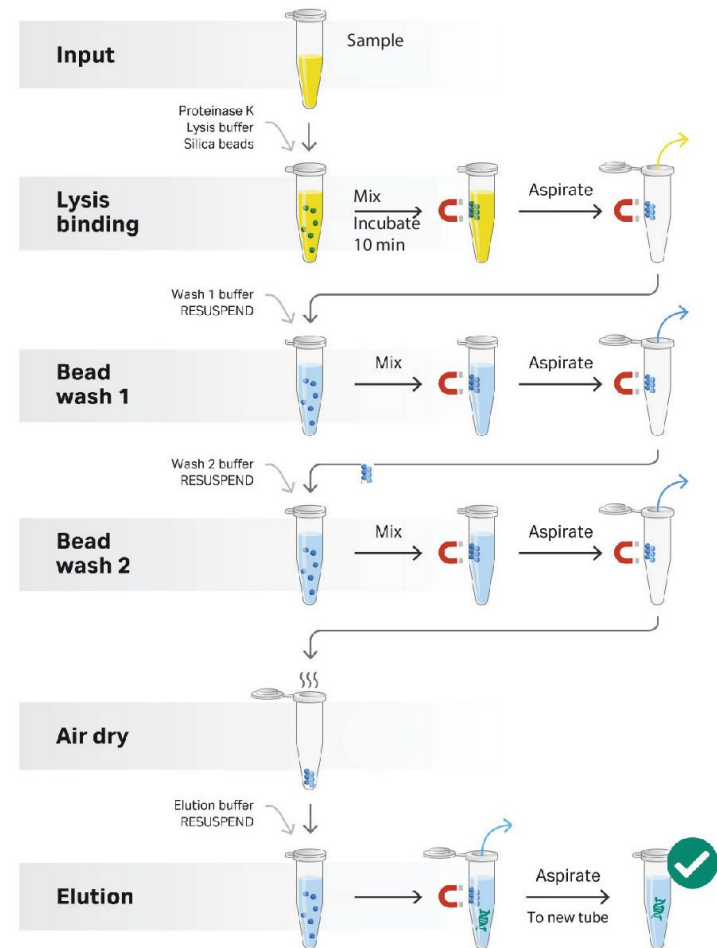


Fig 1. Sera-Xtracta Virus/Pathogen kit workflow

Biomek i7 Dual Hybrid Workstation Automated Cytiva™ Sera-Xtracta Virus/Pathogen Kit

Automated Method Description:

The Cytiva™ Sera-Xtracta Virus/Pathogen kit on Biomek i7 Dual-Hybrid automated workstation method allows the user to isolate viral/bacterial nucleic acids from up to 96 swab samples at a time. The method uses multichannel plate stamping to avoid column effects. It requires only one user interaction should users incubate their samples via the integrated Inheco Deep-Well Incubator or on-deck Inheco Thermoshake AC Peltier.

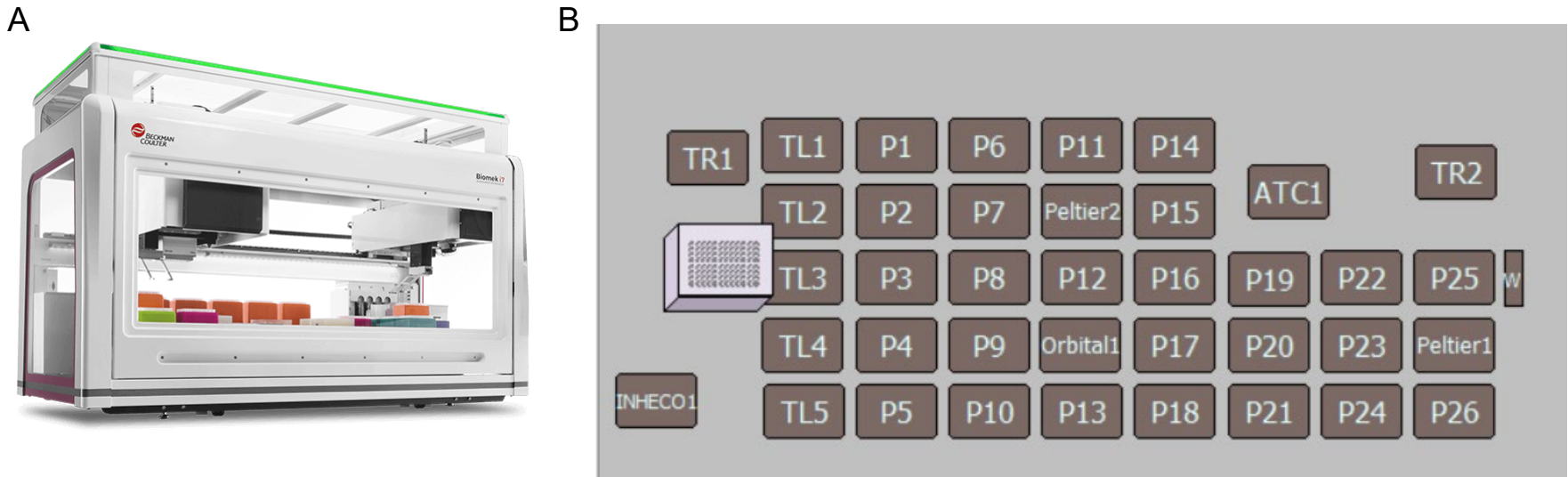


Fig 2.. A. Biomek i7 Dual-Hybrid automated workstation. B. Diagram of the deck layout employed in this method.

Biomek i7 Dual Hybrid Workstation Automated Cytiva™ Sera-Xtracta Virus/Pathogen Kit

Major Process	24 Samples	48 Samples	96 Samples
Instrument Setup	30 min	30 min	30 min
Reagent Aliquoting	13 min, 30 Sec	24 min, 38 sec	34 min, 21 sec
Lysis Binding	21 min, 24 sec	21 min, 24 sec	21 min, 24 sec
Wash 1	16 min, 3 sec	16 min, 3 sec	16 min, 3 sec
Wash 2	13 min, 49 sec	13 min, 49 sec	13 min, 49 sec
Elution	5 min, 41 sec	5 min, 41 sec	5 min, 41 sec
Total	1hr, 40 min, 27 sec	1 hr, 52 min, 35 sec	2hr, 1min, 18 sec

Table 1. Time of completion for 24, 48, and 96 sample inputs.

Biomek i7 Dual Hybrid Workstation Automated Cytiva™ Sera-Xtracta Virus/Pathogen Kit

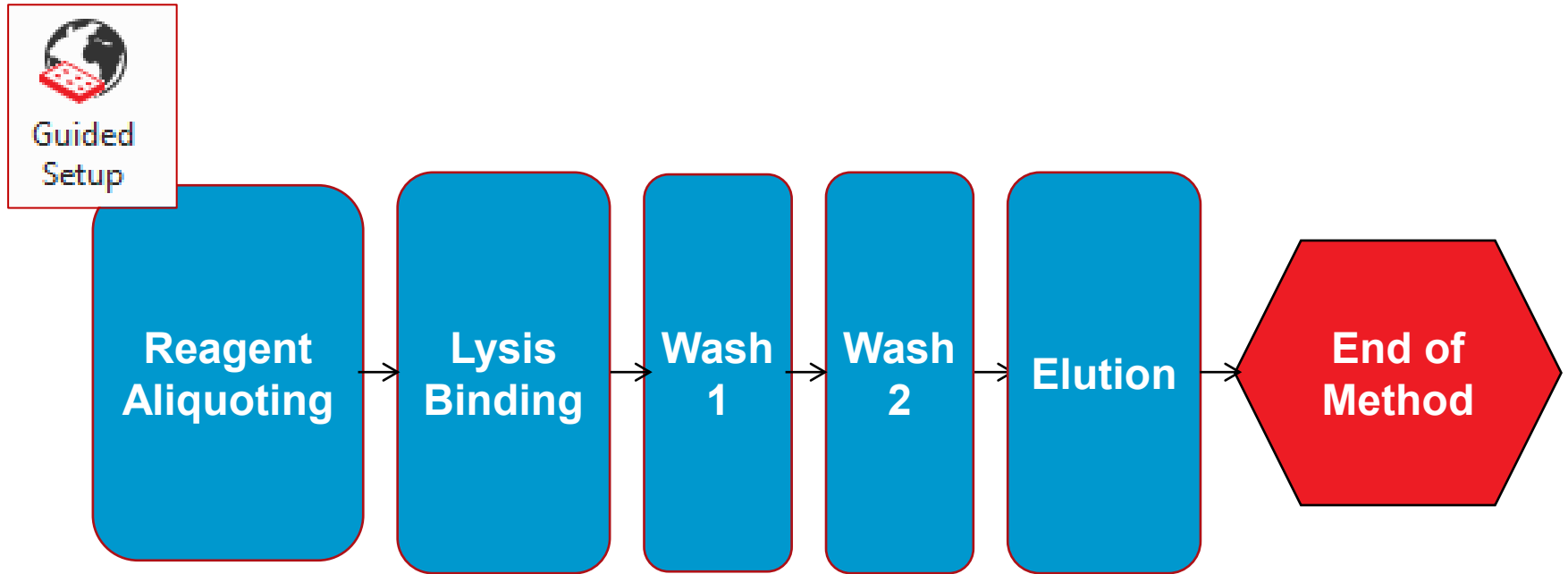


Fig 3. Sera-Xtracta Virus/Pathogen automated workflow.

Biomek i7 Dual Hybrid Workstation Automated Cytiva™ Sera-Xtracta Virus/Pathogen Kit

Demonstrated Method Interface Makes it Easy to Run

Biomek Method Launcher

1. Makes method available outside of the editor
2. Protects against accidental changes from multiple users

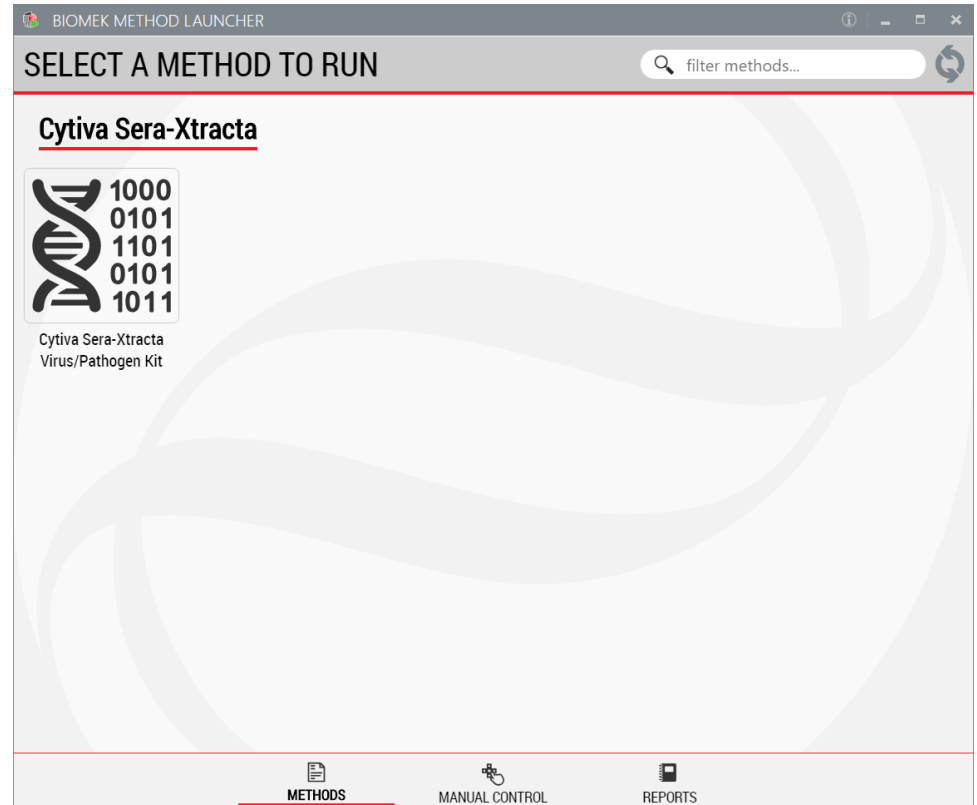


Fig 4. Method selection in the Biomek Method Launcher.

Biomek i7 Dual Hybrid Workstation Automated Cytiva™ Sera-Xtracta Virus/Pathogen Kit

Demonstrated Method Interface Makes it Easy to Run

Method Option Selector (MOS)

Dynamic, HTML-driven interface allows the user to select a number of options to configure the workflow as desired.

The screenshot displays the Method Option Selector (MOS) interface for the Beckman Coulter Cytiva Sera-Xtracta Virus/Pathogen Kit. The interface is titled "Beckman Coulter® Cytiva™ Sera-Xtracta Virus/Pathogen Kit" and is optimized for Biomek i-Series workstations, automated by Beckman Coulter, Inc. The "Method Options" section includes three input fields: "Enter Number of Samples" with a value of 96 (range 8-96 samples), "Enter the Sample Volume in Microliters" with a value of 400 (range 100-400 microliters), and "Incubation Method" with a dropdown menu set to "Incubate on-deck with Inheco Incubator". At the bottom, there are two buttons: "Start run" (green play icon) and "Abort" (red stop icon). The Beckman Coulter logo is visible in the bottom right corner.

Fig 5. Method Option Selector for the Cytiva Sera-Xtracta Virus/Pathogen kit automated method.

Biomek i7 Dual Hybrid Workstation Automated Cytiva™ Sera-Xtracta Virus/Pathogen Kit

Demonstrated Method Interface Makes it Easy to Run

MOS options that enable flexibility, reduce errors and provide greater walk-away time

- 1. Enter Number of Samples:** The Cytiva™ Sera-Xtracta Virus/Pathogen automated method accepts 8-96 samples in multiples of 8 (one column of the plate).
- 2. Specify the sample input volume:** Set the samples' volumes from 100µL to 400µL. The sample volume in each well must be uniform.
- 3. Select the desired process for incubating the sample plate during lysis:** Choose from three different incubation options:
 - 1. Incubate on-deck:* Integrated Inheco Deep Well Single Plate Incubator
 - 2. Incubate on-deck:* An on-deck Inheco Thermoshake AC Peltier
 - 3. Incubate off-deck.* Pause the system at the incubation step to allow the user to remove the plate from the deck and manually incubate. The user replaces the plate to the deck after incubation is complete.

Biomek i7 Dual Hybrid Workstation Automated Cytiva™ Sera-Xtracta Virus/Pathogen Kit

Demonstrated Method Interface Makes it Easy to Run

Guided Labware Setup and Reagent Calculation

- Provides clear instructions for setup
- Reduces setup errors
- GLS diagrams may be printed or saved as a PDF for future reference.

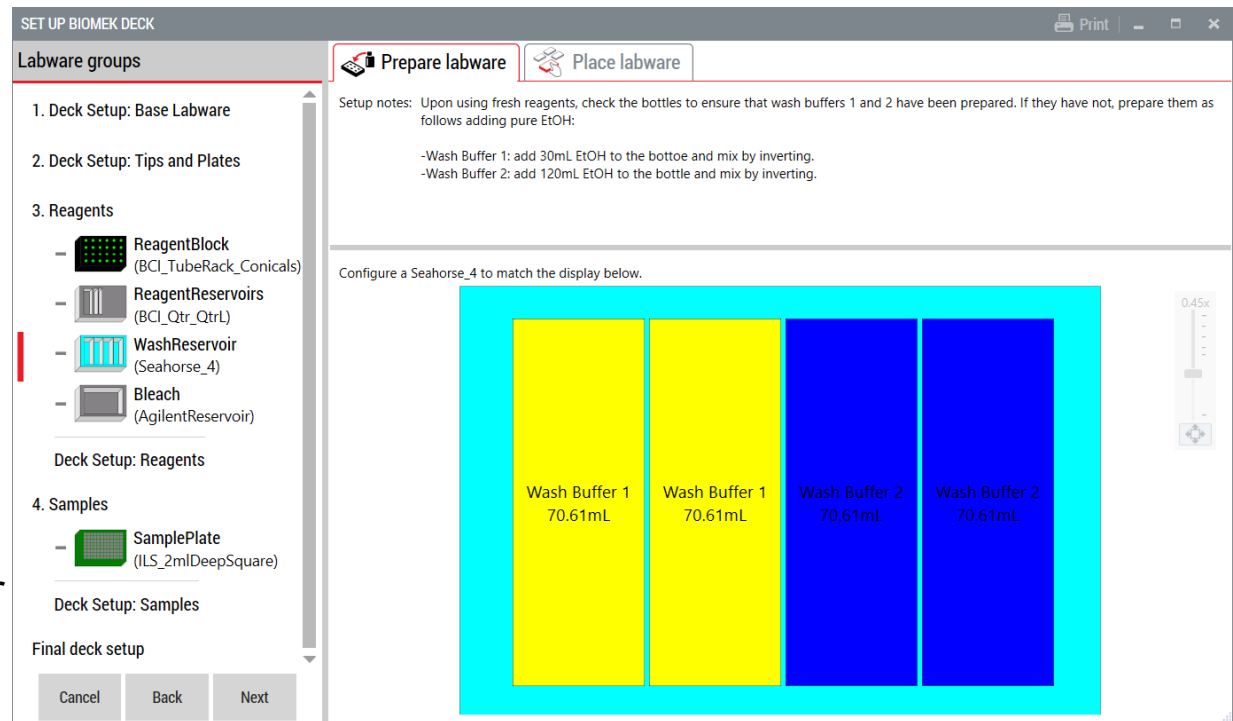


Fig 6. Example of visualized reagent preparation instructions within the Guided Labware Setup steps.

Biomek i7 Dual Hybrid Workstation Automated Cytiva™ Sera-Xtracta Virus/Pathogen Kit

Experimental Details:

All reagent volumes and preparations were prepared according to the method IFU. Samples of HEK 293 cells in a concentration of 3×10^5 /mL in 200µL of viral transport media were spiked with heat-inactivated SARS-CoV-2 RNA for a final concentration of 1 copy/µL. The samples were then processed with the same workflow but different incubation methods for 10 min at 60°C for lysate binding; some samples (n=10) were incubated via the integrated Inheco Deep Well Incubator, while other samples (n=15) were incubated via the on-deck Inheco Thermoshake AC Peltier.

Following elution of the extracted nucleic acid the ct scores of each sample were measured by a QuantStudio™ 6 Flex. The samples were prepared with TaqMan Reagents, Bio-Rad 4x Reliance One-Step Master Mix, and IDT 2019-nCoV_N1 primer.

Biomek i7 Dual Hybrid Workstation Automated Cytiva™ Sera-Xtracta Virus/Pathogen Kit

Samples processed by either the Inheco incubator or the on-deck Peltier both provided Ct means below 40 with minimal deviation.

Samples	Ct mean	St.Dev	Incubation Method
15	35.961	0.652	Peltier
10	36.264	0.462	Inheco Incubator

Table 2. Ct mean and standard deviations and number of samples tested for both incubation methods.

Biomek i7 Dual Hybrid Workstation Automated Cytiva™ Sera-Xtracta Virus/Pathogen Kit

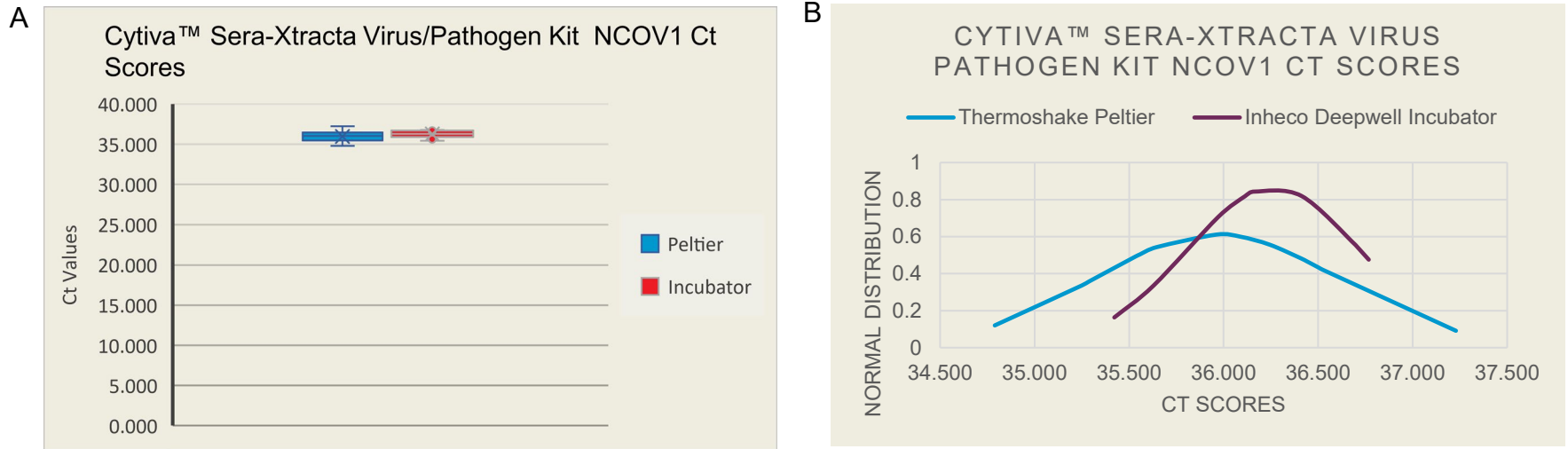


Fig 7. Measured Ct values by incubation method. A. Ct score ranges. B. Ct score normal distributions.

Intellectual Property

This document contains confidential and proprietary information owned by Beckman Coulter, Inc.

Any disclosure, use, copying or distribution of this information without the express written consent of Beckman Coulter, Inc. is strictly prohibited.