



Biomek i5 Nucleic Acid Extraction Solution – Viral Nucleic Acid

RNAadvance Viral and RNAadvance Viral XP



Methods

The method performs extraction of nucleic acid from nasopharyngeal (NP) or oropharyngeal (OP) swabs in transport media, and saliva in a 96-well plate format. Starting input volumes of 200 μ L from a sample are lysed followed by nucleic acid paramagnetic bead binding. The bound nucleic acid and paramagnetic beads isolated are then captured with a magnet and the residual solution is transferred to waste. The beads are washed to remove contaminants and the viral nucleic acids are eluted from the beads in nuclease-free water.

Extraction Process Workflow

RNAadvance Viral Workflow



RNAadvance Viral XP Workflow



Total Estimated Time

		RNAadvance Viral	RNAadvance Viral XP
96 samples (1 plate)	Hands-on Time	15 min	10 min
	Total Time	1 hr, 15 min	55 min
192 samples (2 plates)	Hands-on Time	15 min	10 min
	Total Time	1 hr, 30 min	1 hr, 15 min

Input Material, Reagents, Consumables

Input Material

Sample Type	Transport Media	RNAadvance Viral	RNAadvance Viral XP	Storage Temperature
NP/OP swabs	VTM/UTM	200 μ L	200 μ L	2 - 8 °C < 72 hrs or -70 °C > 72hrs
		200 μ L	NR	Outside collection device recommend storage condition
Saliva	none	200 μ L	NR	Ambient < 72 hr

NR = Not Recommended

Reagents

Description	Supplier	Storage Temperature
RNAAdvance Viral Reagent Kit, 768 Preps	Beckman Coulter	Proteinase K -15 to -25°C Other Components 15 – 30°C
RNAAdvance Viral XP Reagent Kit, 1056 Preps	Beckman Coulter	15 – 30°C
100% Ethanol (Molecular Grade)	User	Room Temp
100% Isopropanol (Molecular Grade)	User	Room Temp
Nuclease-free water (Molecular Grade)	User	Room Temp

Consumable Plastics

Description	Supplier	RNAAdvance Viral Quantity - 2 plates	RNAAdvance Viral XP Quantity - 2 plates
1025µL Pipette Tips, non-Sterile Filtered	Beckman	6	-
190 µL Pipette Tips, non-Sterile Filtered	Beckman	3	6
Quarter Reservoir	Beckman	1	-
Quarter Reservoir, Divided by Length	Beckman	1	-
Reservoir, Half	Beckman	1	-
Hard-Shell® Thin-Wall 96-well Skirted PCR Plates, clear wells	BioRad	3	2
V-bottom 2.2 mL plate. 96-well. PP Sterile	ABgene	7	4
Reservoir 96 Well Pyramid PP 287 mL 25/CS	Agilent	3	3

Method Overview

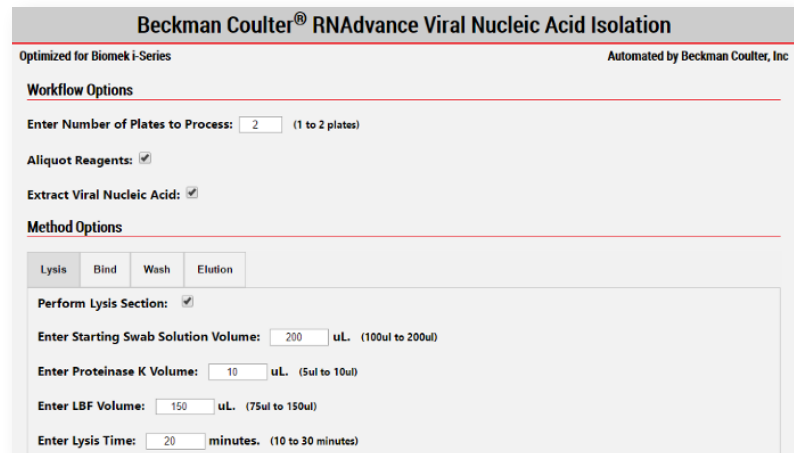
Biomek Method Launcher (BML)

BML organizes methods into useful groups.



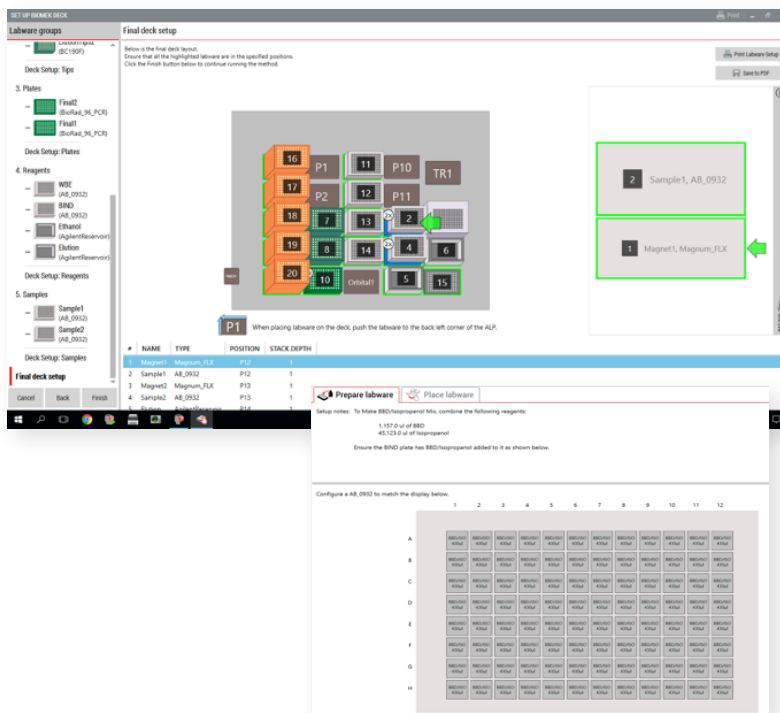
Method Options Selector (MOS)

MOS enables flexibility specific to your sample process batch size, process options and workflow customization.



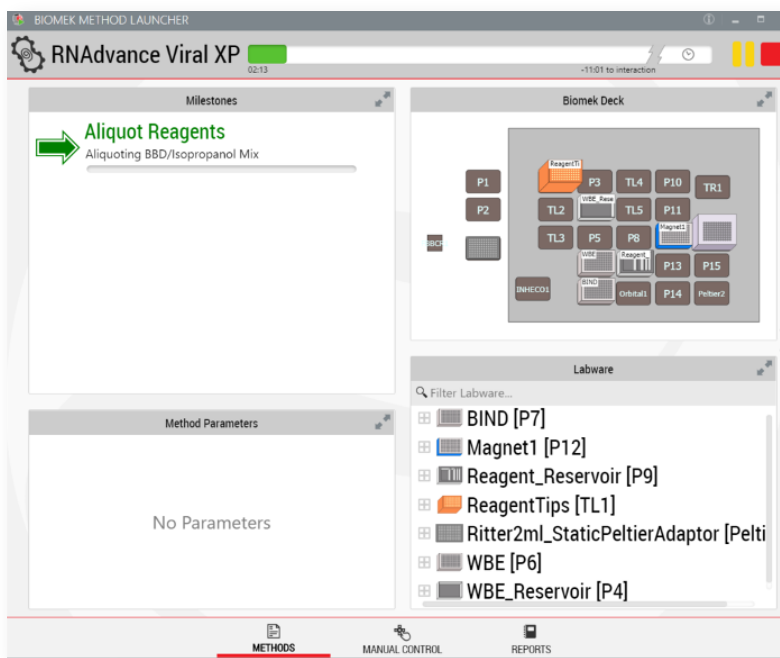
Guided Labware Setup (GLS)

GLS provides clear instructions to set up the instrument deck with calculated reagent volume and step-by-step instructions to prepare reagents based on options selected.



Run Status Screen (RSS)

RSS shows the progress of run, current activity and time to completion.



For Research Use Only. Not intended or validated for use in the diagnosis of disease or other conditions.

©2020 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 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. Not intended or validated for use in the diagnosis of disease or other conditions. This protocol is for demonstration only, and is not validated by Beckman Coulter.



For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at [beckman.com](https://www.beckman.com)

AAG-742707.20