

Purification of Viral RNA from Sputum with the Maxwell[®] HT Viral TNA Kit, Custom

Purify viral RNA from sputum using the Maxwell[®] HT Viral TNA Kit, Custom with the KingFisher[™] Flex Purification System.

Kit:	Maxwell [®] HT Viral TNA Kit, Custom (Cat.# AX2340)	This protocol was developed by
Analyses:	RT-qPCR for detection of Respiratory Syncytial Virus (RSV) and Influenza B.	Promega Applications Scientists and is intended for research use only. Users are responsible for
Sample Type(s):	Sputum	determining suitability of the protocol for their application.
Input:	200µl	For further information, contact Technical Services at: techserv@promega.com
Materials Required:		

- DTT, Molecular Grade (Cat.# V3151)
- Nuclease-Free Water (Cat.# P1193)
- PBS, pH 7.2 (Gibco Cat.# 20012027) or similar
- Maxwell[®] HT Viral TNA Kit, Custom (Cat.# AX2340)
- 4/40 Wash Solution (Cat.# A2221)
- Alcohol Wash, Blood (Cat.# MD1411)
- 80% Ethanol
- 100% Isopropanol
- KingFisher[™] Flex Purification System (ThermoFisher Scientific, Cat.# 24074431)
- KingFisher Deep-well 96 Plate (ThermoFisher Scientific, Cat.# 95040450)
- KingFisher 96 tip comb for DW magnets (ThermoFisher Scientific, Cat.# 97002534)
- KingFisher[™] Flex Run Protocol (Maxwell_HT_Viral_TNAv1_2_RT Elution.bdz)

Sputum Processing¹:

- 1. Weigh appropriate amount of DTT, and rehydrate in Nuclease-Free Water to 500mM final concentration. Mix gently by pipetting to dissolve. DTT must be freshly made.
- 2. Prepare a 1:51 dilution of the 500mM DTT in PBS, pH 7.2. For example, add 100μl of 500mM DTT to 5.0ml of PBS.
- 3. Add an equal volume of diluted DTT in PBS to the sputum sample.
- 4. Incubate at room temperature with intermittent mixing by inversion until liquified or for a maximum of 30 minutes.
- 5. Proceed with purification.



Purification Protocol:

- 1. Prepare 4/40 Wash Buffer and Alcohol Wash Buffer as indicated on the bottles.
- 2. Prepare KingFisher[™] plates:
 - a. Tip Plate: Add Tip comb for KingFisher Deep-Well plate.
 - b. Elution Plate: Add 110µl of Nuclease-Free Water per well.
 - c. 4_40 Wash 1: Add 900µl of 4/40 Wash Buffer per well.
 - d. Alcohol Wash 2: Add 450µl of Alcohol Wash Buffer per well.
 - e. Ethanol Wash 3: Add 450µl of 80% ethanol per well.
 - f. Lysis and Bind: Add the following reagents to each well*:
 - i. 200µl of Lysis Buffer
 - ii. 35µl of Proteinase K
 *Optional: A master mix of the above reagents (plus at least 10% overage) can be prepared immediately before sample addition and 235µl added to each well.
- 3. Transfer 200µl of DTT-treated sputum to each well of the Lysis and Bind Plate.
- 4. Start the KingFisher[™] Flex Run Protocol (Maxwell_HT_Viral_TNAv1_2_RT Elution.bdz).
- 5. Load the KingFisher[™] 96 Deep Well Plates as directed by the instrument software.
- 6. After the heated lysis step, add 530µl of 100% Isopropanol and 35µl of MagneSil[®] RED resin (vortex vigorously to resuspend prior to addition)[^].
 ^Optional: A master mix of the above reagents (plus at least 10% overage) can be prepared and 565µl added to each well. Vortex master mix vigorously before adding to wells.
- 7. Continue the KingFisher[™] Flex Run Protocol until complete.



Product Application



Results:

Detection of RSV and Influenza B RNA extracted from sputum. Sputum was treated with diluted DTT in PBS for 30 minutes. RSV A and Influenza B (Hong Kong) virus were reconstituted from Helix Elite™ Inactivated Standard Inactivated Influenza A/B and Respiratory Syncytial Virus (Microbiologics Cat.# HE0044N) and spiked into treated sputum. High virus sample contains approximately 2 x 10⁵ copies of Influenza B and RSV A per 200µl sample. Low virus sample is a 1:100 dilution of the high virus sample in treated sputum. 200µl of the spiked sputum was processed with the Maxwell® HT Viral TNA Kit, Custom on the KingFisher™ Flex Purification System as described above. Following nucleic acid purification, presence of RSV A and Influenza B was detected by RT-qPCR using GoTaq® 1-Step Probe qPCR System (Cat.# A6121). Each reaction contained 5µl of eluate with 12.5µl of the GoTaq® Probe qPCR Master Mix with dUTP, 0.5µl of GoScript™ RT Mix for 1-Step RT-qPCR, 1000nM forward and reverse primers and 200nM probe for RSV² or Influenza B³, and Nuclease-Free Water added to a final volume of 25µl. 1-step RT-qPCR thermal cycling was as follows³: reverse transcription at 50°C for 30 minutes, hot-start activation at 95°C for 2 minutes, and then 45 cycles of denaturation at 95°C for 15 seconds and annealing/extension at 55°C for 30 seconds, with signal acquisition during the annealing/extension stage of cycling. Data represent the average of duplicate purifications amplified in duplicate. Error bars indicate the standard deviation. Influenza B was not detected in the RNA extracted from the low virus sample.

References:

- Processing of Sputum Specimens for Nucleic Acid Extraction, Centers for Disease Control <u>https://www.cdc.gov/coronavirus/2019-ncov/downloads/processing-sputum-specimens.pdf</u> Accessed 3/12/2020.
- 2. Fry, A.M., *et al.*, (2010) The Burden of Hospitalized Lower Respiratory Tract Infection due to Respiratory Syncytial Virus in Rural Thailand, *PLoS One*. *5*, e15098.
- 3. Selvaraju, S.B., *et al.*, (2010) Evaluation of Three Influenza A and B Real-Time Reverse Transcription-PCR Assays and a New 2009 H1N1 Assay for Detection of Influenza Viruses, *Journal of Clinical Microbiology.* **48**, 3870-3875.