

Automated Zika Virus RNA Isolation from Biological Liquids.

Isolate Zika Virus RNA from plasma and other biological liquids using the Maxwell® RSC Instrument.

Kit: Maxwell® RSC Viral Total Nucleic Acid Purification Kit (Cat. #AS1330)

Analyses: RT-qPCR

Sample Type(s): Plasma, Blood, Saliva

Input: up to 300µl of Sample

Materials Required:

- Maxwell® RSC Total Nucleic Acid Purification Kit (Cat. #AS1330)
- Maxwell® RSC Instrument (Cat. #AS4500)
- Heat Block capable of 56°C
- Microcentrifuge tubes

This protocol was developed by Promega Applications Scientists and is intended for research use only.

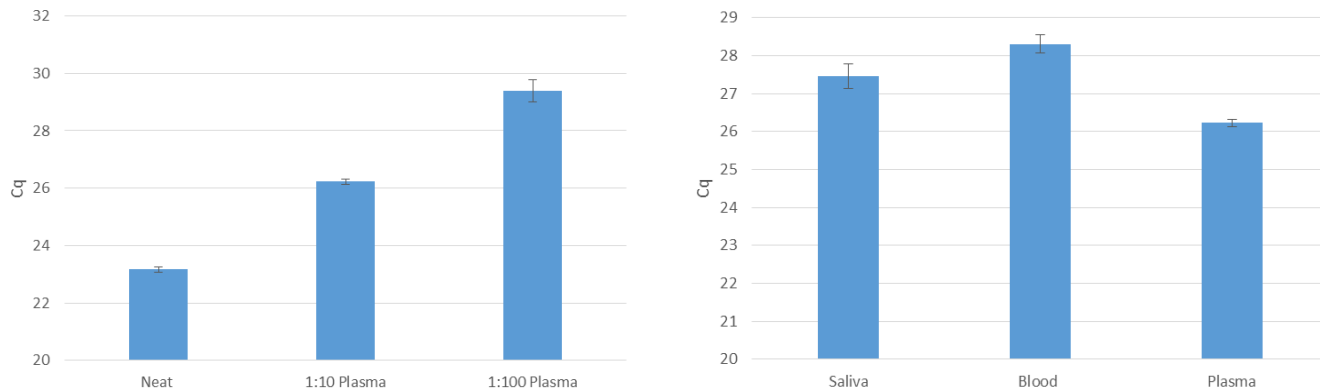
Users are responsible for determining suitability of the protocol for their application.

Further information can be found in Technical Manual #TM420, available at: www.promega.com/protocols or for further information, please contact techserv@promega.com

Protocol:

1. Add 20µl Proteinase K and 200µl Lysis Buffer to up to 300µl of sample.
2. Vortex for 10 seconds.
2. Incubate samples at 56°C for 10 minutes at 56°C.
3. Load sample into well #1 of the Maxwell Cartridge and place a Maxwell plunger into well #8. Add 50µl of Nuclease-Free Water to the elution tube.
4. Select the RSC Viral TNA method for processing the samples on the Maxwell® RSC Instrument.

Example Performance of Maxwell RSC Total Nucleic Acid Purification Kit with Zika Virus



C_q values from RT-qPCR for Zika Virus with RNA from Plasma (Left). Zika Virus obtained from Zeptomatrix was used for testing (Zeptomatrix, Cat. # NATZIKV-ERCM). Zika Virus was spiked into human plasma at a ratio of 1:10 or 1:100 v/v. Viral recovery was tested by amplification with GoTaq® 1-step RT-qPCR System using Zika virus specific primers. n=3 for each condition.

C_q values from RT-qPCR for Zika Virus with RNA from Plasma, Blood, and Saliva (Left). Zika Virus obtained from Zeptomatrix was used for testing (Zeptomatrix, Cat. # NATZIKV-ERCM). Zika Virus stock was spiked into whole blood, plasma, or saliva at 1:10 v/v ratio. Viral recovery was tested by amplification with GoTaq® 1-step RT-qPCR System using Zika virus specific primers. n=3 for each condition.