

# MagMAX Viral/Pathogen kit protocol changes to enable increased SARS-CoV-2 testing throughput

## Reduction in sample input volume and wash steps, plus alternative KingFisher plastics

This application note provides performance data on the Applied Biosystems™ MagMAX™ Viral/Pathogen Nucleic Acid Isolation Kit and MagMAX™ Viral/Pathogen II Nucleic Acid Isolation Kit, and guidance on the following technical elements for SARS-CoV-2 testing:

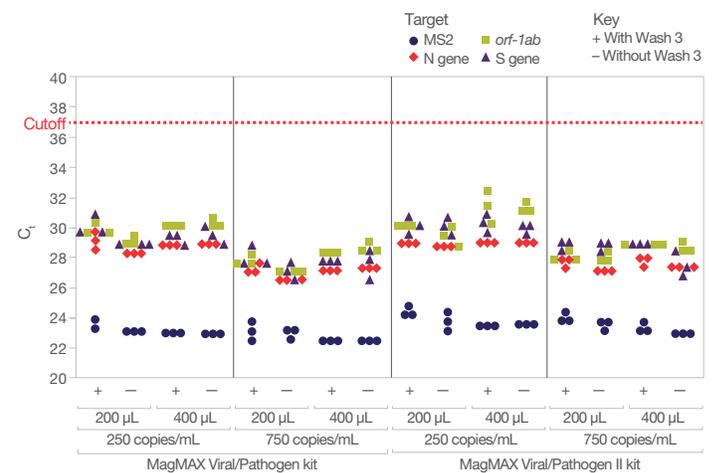
- Reduction of sample input volume from 400  $\mu\text{L}$  to 200  $\mu\text{L}$
- Reduction from 3 wash steps to 2 wash steps
- Use of alternative plastic plates and tip combs for Thermo Scientific™ KingFisher™ instruments

### Equivalent performance of MagMAX Viral/Pathogen and MagMAX Viral/Pathogen II Nucleic Acid Isolation Kits

The MagMAX Viral/Pathogen kit (Cat. No. A48310) and MagMAX Viral/Pathogen II kit (Cat. No. A48383R) demonstrate equivalent performance as measured by  $C_t$  values of spiked SARS-CoV-2 RNA, enabling greater flexibility and testing capacity (Figure 1). The protocols for the MagMAX Viral/Pathogen kit and MagMAX Viral/Pathogen II kit result in equivalent RNA purification efficiency whether using 400  $\mu\text{L}$  or 200  $\mu\text{L}$  of sample input (and normalized going into the qRT-PCR reaction). The component volumes per well and  $C_t$  values are shown in Table 1 and Figure 1, respectively. Samples were processed on a Thermo Scientific™ KingFisher™ Flex Purification System with 96 Deep-Well Head (Cat. No. 5400630).

**Table 1. Sample prep protocol for MagMAX Viral/Pathogen kits.**

	400 $\mu\text{L}$ sample input	200 $\mu\text{L}$ sample input
Component	Volume per well	
Proteinase K	10 $\mu\text{L}$	5 $\mu\text{L}$
Sample (or nuclease-free water for negative control)	400 $\mu\text{L}$	200 $\mu\text{L}$
Binding bead mix (binding solution + beads)	550 $\mu\text{L}$ (530 $\mu\text{L}$ + 20 $\mu\text{L}$ )	275 $\mu\text{L}$ (265 $\mu\text{L}$ + 10 $\mu\text{L}$ )
MS2 phage control	10 $\mu\text{L}$	5 $\mu\text{L}$



**Figure 1.  $C_t$  values for contrived SARS-CoV-2-positive nasopharyngeal swab samples with different sample preparation workflows.** Nasopharyngeal swab specimens were spiked with 250 copies/mL or 750 copies/mL of SARS-CoV-2 viral RNA (1x or 3x the limit of detection, respectively), and RNA was extracted from 200  $\mu\text{L}$  or 400  $\mu\text{L}$  of the spiked sample in triplicate using either the MagMAX Viral/Pathogen or MagMAX Viral/Pathogen II kit with or without Wash 3. After elution in 50  $\mu\text{L}$ , the samples were assayed by qRT-PCR. All negative samples had  $C_t$  values of “Undetected” (data not shown).

### Reduction from 3 wash steps to 2 wash steps

The MagMAX Viral/Pathogen kit and MagMAX Viral/Pathogen II kit protocols for SARS-CoV-2 testing using the KingFisher Flex Purification System with 96 Deep-Well Head support both 3 wash steps (Tables 2 and 3) and 2 wash steps (Tables 4 and 5).

**Table 2. Protocol for 400 µL sample input with 3 wash steps (script name: MVP\_Flex\_400ul.bdz).**

Plate ID	Plate position	Plate type	Reagent	Volume per well
Wash 1 plate	2	KingFisher deep-well 96 plate	Wash buffer	1,000 µL
Wash 2 plate	3		80% ethanol	1,000 µL
Wash 3 plate	4		80% ethanol	500 µL
Elution plate	5		Elution solution	50 µL
Tip comb	6		Place a KingFisher 96 tip comb for deep-well magnets in a KingFisher 96 microplate or alternative tip comb-presenting plate	

**Table 3. Protocol for 200 µL sample input with 3 wash steps (script name: MVP\_Flex\_200ul.bdz).**

Plate ID	Plate position	Plate type	Reagent	Volume per well
Wash 1 plate	2	KingFisher deep-well 96 plate	Wash buffer	500 µL
Wash 2 plate	3		80% ethanol	500 µL
Wash 3 plate	4		80% ethanol	250 µL
Elution plate	5		Elution solution	50 µL
Tip comb	6		Place a KingFisher 96 tip comb for deep-well magnets in a KingFisher 96 microplate or alternative tip comb-presenting plate	

Updated scripts may be obtained at [thermofisher.com/magmaxviralpathogen](http://thermofisher.com/magmaxviralpathogen), or by editing the existing scripts according to Tables 4 and 5.

**Table 4. Protocol for 400 µL sample input with 2 wash steps (script name: MVP\_2Wash\_400\_Flex.bdz).**

Plate ID	Plate position	Plate type	Reagent	Volume per well	
Wash 1 plate	2	KingFisher deep-well 96 plate	Wash Buffer	1,000 µL	
Wash 2 plate	3		80% ethanol	1,000 µL	
Elution plate	4		Elution solution	50 µL	
Tip comb	5		Place a KingFisher 96 tip comb for deep-well magnets in a KingFisher 96 microplate or alternative tip comb-presenting plate		

**Table 5. Protocol for 200 µL sample input with 2 wash steps (script name: MVP\_2Wash\_200\_Flex.bdz).**

Plate ID	Plate position	Plate type	Reagent	Volume per well	
Wash 1 plate	2	KingFisher deep-well 96 plate	Wash buffer	500 µL	
Wash 2 plate	3		80% ethanol	1,000 µL	
Elution plate	4		Elution solution	50 µL	
Tip comb	5		Place a KingFisher 96 tip comb for deep-well magnets in a KingFisher 96 microplate or alternative tip comb-presenting plate		

Here we show the reduction from 3 wash steps to 2 wash steps without any loss of sensitivity for the MagMAX Viral/Pathogen kit and the MagMAX Viral/Pathogen II kit (Figure 1). This 2-wash protocol conserves both processing plastics and wash reagents as well as time and effort. This protocol has been tested with qRT-PCR; other applications have not been evaluated.

## Alternative plastics for KingFisher instruments

The following products can be used as alternatives for the existing Thermo Scientific™ KingFisher™ 96 deep-well, V-bottom plate (Cat. No. 95040450).

Plate description	KingFisher instrument compatibility	Quantity	Cat. No.
<b>New</b> KingFisher 96 deep-well plate	Flex, Duo Prime	50 plates per case	95040455
<b>New</b> KingFisher 96 deep-well plate			A48424*
<b>New</b> KingFisher 96 deep-well plate	Flex, Duo Prime, Presto		A48305

The following products can be used as alternatives for the existing Thermo Scientific™ KingFisher™ 96 tip comb for deep-well magnets (Cat. No. 97002534).

Tip comb description	KingFisher instrument compatibility	Quantity	Cat. No.
<b>New</b> KingFisher 96 tip comb for deep-well plates	Flex, Presto	100 tip combs per case	A48438
<b>New</b> KingFisher 96 tip comb for deep-well plates		50 tip combs per case	A48414*

## Ordering information

Product	Quantity	Cat. No.
<b>Recommended</b> MagMAX Viral/Pathogen II Nucleic Acid Isolation Kit	Up to 2,000 preps	A48383R
MagMAX Viral/Pathogen Nucleic Acid Isolation Kit	Up to 2,000 preps	A48310
KingFisher Flex Purification System with 96 Deep-Well Head	1 system	5400630
KingFisher 96 deep-well plate, V-bottom, polypropylene	50/box	95040450
<b>New</b> KingFisher 96 deep-well plate, V-bottom, polypropylene	50/box	A48305
<b>New</b> KingFisher 96 deep-well plate (for Flex and Duo Prime instruments only)	50/box	95040455
<b>New</b> KingFisher 96 deep-well plate (for Flex and Duo Prime instruments only)	50/box	A48424*
<b>New</b> KingFisher 96 tip comb for deep-well plates	100/box	A48438
<b>New</b> KingFisher 96 tip comb for deep-well plates	50/box	A48414*
KingFisher 96 (200 µL) microplate	48/box	97002540
KingFisher 96 tip comb for deep-well magnets	100/box	97002534

\* Not available in all regions.

Find out more about MagMAX Viral/Pathogen Nucleic Acid Isolation Kits at [thermofisher.com/magmaxviralpathogen](https://thermofisher.com/magmaxviralpathogen)

Find out more about KingFisher plastics at [thermofisher.com/kingfisherplastics](https://thermofisher.com/kingfisherplastics)

**ThermoFisher**  
SCIENTIFIC