# MagMAX<sup>™</sup> Saliva gDNA Isolation Kit

High throughput isolation of gDNA from saliva

Catalog Numbers A39059, A39060

Pub. No. MAN0017722 Rev. A.0

Â

**WARNING!** Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from **thermofisher.com/support**.

## **Product description**

The Applied Biosystems<sup>™</sup> MagMAX<sup>™</sup> Saliva gDNA Isolation Kit is developed for scalable, rapid purification of high-quality DNA from saliva, both preserved (saliva within a stabilizing reagent or a preservative) and fresh. You can use the DNA purified with this kit in a broad range of molecular biology downstream applications, such as sequencing, genotyping, and qPCR. This protocol guides through automated isolation of DNA from saliva using the KingFisher<sup>™</sup> Flex and the KingFisher<sup>™</sup> Duo Prime.

## **Contents and storage**

Reagents that are provided in the kit are sufficient for 100 reactions (Cat. No. A39059) or 500 reactions (Cat. No. A39060) using small volume ( $\leq$ 500 µL) inputs.

Table 1 Components of MagMAX<sup>™</sup> Saliva gDNA Isolation Kit

Component	100 reactions	500 reactions	Storage
Lysis/Binding Solution	55 mL	275 mL	
gDNA Binding Beads	4.5 mL	22 mL	15°C to 30°C
Wash I Solution	110 mL	2 × 275 mL	
Elution Solution	12 mL	60 mL	

For 1,000 reaction volume, use Cat. No. A39063 (Lysis/Binding Solution), A39061 (gDNA Binding Beads), A39062 (Wash I Solution), and A39064 (Elution Solution).

## Required materials not supplied

Unless otherwise indicated, all materials are available through **thermofisher.com**. MLS: Fisher Scientific (**fisherscientific.com**) or other major laboratory supplier.

Item	Source	
Magnetic particle processor (one of the following, depending on quantity/volume of sample to be processed):		
KingFisher™ Duo Prime Magnetic Particle Processor	5400110	
<i>For small volume sample<sup>[1]</sup>:</i> KingFisher™ Flex Magnetic Particle Processor with 96 Deep-Well Head	5400630	
<i>For large volume sample</i> <sup>[2]</sup> : KingFisher <sup>™</sup> Flex Magnetic Particle Processor with 24 Deep-Well Head	5400640	
Equipment		
Adjustable micropipettors	MLS	
Multi-channel micropipettors	MLS	
Consumables		
Deep-well plates:		
<i>For small volume sample</i> <sup>[1]</sup> : KingFisher™ Deepwell 96 Plate	95040450	
<i>For large volume sample</i> <sup>[2]</sup> : KingFisher <sup>™</sup> Flex 24 Deep- Well Plates	95040470	
96-well standard plates (for use with KingFisher™ Flex only; tip comb placement and elution plate and/or eluate storage):		
KingFisher™ 96 KF microplate	97002540	
Tip comb, compatible with the magnetic particle processor used:		
KingFisher™ Duo Prime 12-tip comb, for use with KingFisher™ deep-well 96 plate	97003500	
KingFisher™ Duo Prime 6 tip comb, for use with the Flex 24 deep-well plate	97003510	
KingFisher <sup>™</sup> 96 tip comb for DW magnets, for Flex protocol only	97002534	
KingFisher™ Flex 24 Deep Well Tip Comb and plate, for Flex protocol only	97002610	
Materials		
MicroAmp <sup>™</sup> Clear Adhesive Film	4306311	

Item	Source
Reagents	
Ethanol, 96–100% (molecular biology grade)	MLS
Nuclease-free Water	AM9932

<sup>[1]</sup> Small volume saliva is ≤500 µL.

<sup>[2]</sup> Large volume saliva is 500 µL-2 mL.

## General guidelines

- Perform all steps at room temperature (20–25°C), unless otherwise noted.
- Precipitates can occur if the Lysis/Binding Solution is stored when room temperature is too cold. If there are precipitates, warm the Lysis/Binding Solution at 37°C and gently mix to dissolve the precipitates. Avoid creating bubbles.
- · Per-plate volumes for reagent mixes are sufficient for one plate plus overage. To calculate volumes for other sample numbers, see the per-well volume and add 10% overage.
- (Optional): To prevent evaporation and contamination, cover the prepared processing plates with paraffin film or MicroAmp<sup>™</sup> Clear Adhesive Film until they are loaded into the instrument.
- When isolating from preserved or stabilized saliva, verify with manufacturer recommendations before processing. Many saliva collection devices require an upfront incubation before sample processing can occur. Failure to do so can affect vields.

## Guidelines for DNA Lysis/Binding Bead Mix

- Vortex gDNA Binding Beads thoroughly, then combine them with Lysis/Binding Solution in a nuclease-free tube and invert the tube until homogeneous. You can store this mixture for up to 1 day before aliquoting into the plates.
- Ensure that the beads stay fully mixed within the solution during pipetting.
- Avoid creating bubbles during mixing and aliquoting.

## Before first use of the kit

Prepare Wash II Solution: Make 80% Ethanol from 100% absolute Ethanol and Nuclease-Free Water.

## Before each use of the kit

Vortex gDNA Binding Beads to fully resuspend the beads before each use.

## Perform DNA purification using KingFisher<sup>™</sup> Flex (small volume: ≤500 µL)

Set up the instrument 1

Ensure that the instrument is set up with the proper magnetic head and the proper heat block, as a. indicated in the following table.

Component	Туре
Magnetic head	96 deep-well magnetic head
Heat block	96 well standard heat block

**IMPORTANT!** Failure to use the proper magnetic head and heat block results in lower yields.

b. Ensure that the proper program (MagMAX\_Saliva\_500µL\_FLEX) has been downloaded from the product page and loaded onto the instrument.

#### Set up the processing **~**

Set up the Wash, Elution, and Tip Comb Plates outside the instrument according to the following table.

occup inc	· P
plates	

Plate ID	Plate position	Plate type	Reagent	Volume per well
≤500 µL saliva input				
Wash I Solution Plate	2	Deep-well	Wash I Solution	1,000 µL
Wash II Solution Plate 1	3	Deep-well	Wash II Solution	1,000 µL
Wash II Solution Plate 2	4	Deep-well	Wash II Solution	500 μL
Elution Plate	5	Standard	Elution Solution	50–100 μL
Tip Comb	6	Place a 96	Deep-well Tip Comb in	a Standard Plate

Note: Load the plates onto the instrument immediately after the Sample Plate has been prepared.

#### 3 Prepare Sample Plate and purify the DNA

**Prepare Sample Plate and a.** Prepare the DNA Lysis/Binding Bead Mix according to the following table.

Component	Volume per well	Volume per plate
≤500 µL saliva input (96 samples	s per plate)	
Lysis/Binding Solution	460 µL	48.58 mL
DNA Binding Beads	40 µL	4.22 mL
Total volume	500 μL	52.80 mL

**b.** Transfer the appropriate amount of the saliva sample to the appropriate wells of a new deep-well plate.

This plate is the Sample Plate.

c. Add 500 µL of the DNA Lysis/Binding Bead Mix to each sample in the Sample Plate.

Note:

- Remix the DNA Lysis/Binding Bead Mix frequently during pipetting to ensure even distribution of beads to all samples or wells. The mixture containing the Binding Beads is viscous. Therefore, pipet slowly to ensure that the correct amount is added.
- After you have added all components, proceed immediately to the instrument processing. There is no need for manual mixing beforehand.

IMPORTANT! Avoid creating bubbles during mixing and aliquoting.

- d. Select the program MagMAX\_Saliva\_500µL\_FLEX on the instrument.
- e. Start the run, then load the prepared plates into position when prompted by the instrument.
- f. At the end of the run, immediately remove the plate from the instrument, then cover the plate or transfer the eluate to the final tube or plate of choice for final storage.

The purified DNA is ready for immediate use. Alternatively, store the plate at -20 °C for long-term storage.

## Perform DNA purification using KingFisher<sup>™</sup> Flex (large volume: 500 µL to 2 mL)

1

Set up the instrument

**a.** Ensure that the instrument is set up with the proper magnetic head and the proper heat block, as indicated in the following table.

Component	Туре
Magnetic head	24 deep-well magnetic head
Heat block	24 well standard heat block

IMPORTANT! Failure to use the proper magnetic head and heat block results in lower yields.

**b.** Ensure that the proper program (**MagMAX\_Saliva\_2mL\_FLEX**) has been downloaded from the product page and loaded onto the instrument.

### 2 Set up the processing plates

Set up the Wash, Elution, and Tip Comb Plates outside the instrument according to the following table and sample input volume.

Plate ID	Plate position	Plate type	Reagent	Volume per well
500 µL – 1 mL saliva inpo	ut			
Wash I Solution Plate 1	2	Deep-well	Wash I Solution	1,000 µL
Wash I Solution Plate 2	3	Deep-well	Wash I Solution	1,000 µL
Wash II Solution Plate 1	4	Deep-well	Wash II Solution	2,000 µL
Wash II Solution Plate 2	5	Deep-well	Wash II Solution	500 μL
Elution Plate	6	Deep-well	Elution Solution	200 µL
Tip Comb	7	KingFisher <sup>™</sup> Flex 24 Deep-well Tip Comb and Plate		
1 mL – 2 mL saliva input				
Wash I Solution Plate 1	2	Deep-well	Wash I Solution	2,000 µL
Wash I Solution Plate 2	3	Deep-well	Wash I Solution	2,000 µL
Wash II Solution Plate 1	4	Deep-well	Wash II Solution	2,000 µL
Wash II Solution Plate 2	5	Deep-well	Wash II Solution	500 μL
Elution Plate	6	Deep-well	Elution Solution	200 µL
Tip Comb	7	KingFisher™ Flex 24 Deep-well Tip Comb and Plate		

Note: Load the plates onto the instrument immediately after the Sample Plate has been prepared.

# **Prepare Sample Plate and** a. Prepare the DNA Lysis/Binding Bead Mix according to the following table and sample input volume.

Component	Volume per well	Volume per plate			
500 μL – 1 mL saliva input (24 sa	500 μL – 1 mL saliva input (24 samples per plate)				
Lysis/Binding Solution	920 μL	24.29 mL			
DNA Binding Beads	80 µL	2.11 mL			
Total volume	1,000 µL	26.40 mL			
1.1 mL – 2 mL saliva input (24 sa	mples per plate)				
Lysis/Binding Solution	1,840 µL	48.58 mL			
DNA Binding Beads	160 µL	4.22 mL			
Total volume	2,000 µL	52.80 mL			

**b.** Transfer the appropriate amount of the saliva sample to the appropriate wells of a new deep-well plate.

This plate is the Sample Plate.

c. Add DNA Lysis/Binding Bead Mix to each sample in the Sample Plate, according to the following table and sample input volume.

For saliva input volume	Add DNA Lysis/Binding Bead Mix
500 μL – 1 mL	1,000 µL
1.1 mL – 2 mL	2,000 µL

2	Prepare Sample Plate
	and purify the DNA
	(continued)

#### Note:

- Remix the DNA Lysis/Binding Bead Mix frequently during pipetting to ensure even distribution of beads to all samples or wells. The mixture containing the Binding Beads is viscous. Therefore, pipet slowly to ensure that the correct amount is added.
- After you have added all components, proceed immediately to the instrument processing. There is no need for manual mixing beforehand.

**IMPORTANT!** Avoid creating bubbles during mixing and aliquoting.

- d. Select the program MagMAX\_Saliva\_2mL\_FLEX on the instrument.
- e. Start the run, then load the prepared plates into position when prompted by the instrument.
- **f.** At the end of the run, immediately remove the plate from the instrument, then cover the plate or transfer the eluate to the final tube or plate of choice for final storage.

The purified DNA is ready for immediate use. Alternatively, store the plate at –20°C for long-term storage.

## Perform DNA purification using KingFisher<sup>™</sup> Duo Prime (small volume: ≤500 µL)

1	Set up the instrument	<b>a</b> . Ensure that the instrument is set up for processing with the proper magnetic head (12 pin) and heat block for your application.			
		<b>b.</b> Ensure that the prop product page and lo	per program ( <b>MagMAX_S</b> aded onto the instrument	aliva_500μL_DUO) has b	een downloaded from the
2	Set up the processing plate	ing Prepare the plate according to the following table. The Sample Row will be prepared in the next section. Table 2 96 Deep-well plate layout (<500 μL saliva input)			
		Row ID	Plate Row	Reagent	Volume per well
		Elution Solution	Α	Elution Solution	50–100 μL
		_	В	Em	npty
		Tip Comb	С	Tip Comb	N/A
		Wash II Solution row 2	D	Wash II Solution	500 μL
		Wash II Solution row 1	E	Wash II Solution	1,000 µL
		Wash I Solution	F	Wash I Solution	1,000 µL
		_	G	Em	npty
		Sample	Н	Sample	Varies

Note: Load the plate onto the instrument immediately after the Sample Row has been prepared.

**3 Prepare Sample Row and a.** Prepare the DNA Lysis/Binding Bead Mix according to the following table.

#### purify the DNA

ComponentVolume per wellVolume per plate<500 μL saliva input (12 samples per plate)</td>Lysis/Binding Solution460 μL6.07 mLDNA Binding Beads40 μL528 μLTotal volume500 μL6.60 mL

**b.** Transfer the appropriate amount of the sample to the appropriate wells (Row H) of the deep-well plate that was previously prepared.

c. Add 500 µL of the DNA Lysis/Binding Bead Mix to each sample in Row H.

#### Note:

- Remix the DNA Lysis/Binding Bead Mix frequently during pipetting to ensure even distribution of beads to all samples or wells. The mixture containing the Binding Beads is viscous. Therefore, pipet slowly to ensure that the correct amount is added.
- After you have added all components, proceed immediately to the instrument processing. There is no need for manual mixing beforehand.

**IMPORTANT!** Avoid creating bubbles during mixing and aliquoting.

d. Select the program MagMAX\_Saliva\_500µL\_DUO on the instrument.

e. Start the run, then load the prepared plate into position when prompted by the instrument.

f. At the end of the run, immediately remove the plate from the instrument, then cover the plate or transfer the eluate to the final tube or plate of choice for final storage.

3	Prepare Sample Row	The purified DNA is ready for immediate use. Alternatively, store the plate at -20°C for long-term
	and purify the DNA	storage.
	(continued)	

## Perform DNA purification using KingFisher<sup>™</sup> Duo Prime (large volume: 500 µL to 2mL)

1	Set up the instrument	<ul><li>a. Ensure that the instruheat block for your a</li><li>b. Ensure that the properties of the product page and load</li></ul>	ument is set up for proces pplication. er program ( <b>MagMAX_S</b> aded onto the instrument	ssing with the proper mag saliva_2mL_DUO) has bee	netic head (6 pin) and en downloaded from the
2	Set up the processing plate	Prepare the 24 deep-well plates according to the following tables and sample input volume. <b>Table 3</b> 24 Deep-well plate layout (500 μL – 1 mL saliva input)			
		Row ID	Plate Row	Reagent	Volume per well
		Plate 1 layout			
		Wash I Solution plate 2	А	Wash I Solution	1,000 µL
		Wash I Solution plate 1	В	Wash I Solution	1,000 µL
		Tip Comb	С	Tip Comb	N/A
		Sample	D	Sample	Varies
		Plate 2 layout			
		Elution Solution	А	Elution Solution	200 µL
		Empty	В		_
		Wash II Solution plate 2	С	Wash II Solution	500 µL
		Wash II Solution plate 1	D	Wash II Solution	2,000 µL
		Table 4 24 Deep-well plate	layout (1 mL – 2 mL saliva	input)	
		Row ID	Plate Row	Reagent	Volume per well
		Plate 1 layout			
		Wash I Solution plate 2	А	Wash I Solution	2,000 µL
		Wash I Solution plate 1	В	Wash I Solution	2,000 µL
		Tip Comb	С	Tip Comb	N/A
		Sample	D	Sample	Varies

**Note:** Load the plates onto the instrument immediately after the Sample Row has been prepared.

А

В

С

D

**Elution Solution** 

Wash II Solution

Wash II Solution

\_

Plate 2 layout Elution Solution

Wash II Solution plate 2

Wash II Solution plate 1

Empty

200 µL

500 µL

2,000 µL

a. Prepare the DNA Lysis/Binding Bead Mix according to the following table and sample input volume.

Component	Volume per well	Volume per plate		
500 μL – 1 mL saliva input (6 samples per plate)				
Lysis/Binding Solution	920 μL	6.07 mL		
DNA Binding Beads	80 µL	528 μL		
Total volume	1,000 µL	6.60 mL		
1.1 mL – 2 mL saliva input (6 samples per plate)				
Lysis/ Binding Solution	1,840 µL	12.14 mL		
DNA Binding Beads	160 μL	1.06 mL		
Total volume	2,000 μL	13.20 mL		

- **b.** Transfer the appropriate amount of the sample to the appropriate wells (Row D of Plate 1) of the deep-well plate that was previously prepared.
- **c.** Add DNA Lysis/Binding Bead Mix to each sample in Row D of Plate 1, according to the following table.

For saliva input volume	Add DNA Lysis/Binding Bead Mix
500 µL – 1 mL	1,000 µL
1.1 mL – 2 mL	2,000 µL

#### Note:

- Remix the DNA Lysis/Binding Bead Mix frequently during pipetting to ensure even distribution of beads to all samples or wells. The mixture containing the Binding Beads is viscous. Therefore, pipet slowly to ensure that the correct amount is added.
- After you have added all components, proceed immediately to the instrument processing. There is no need for manual mixing beforehand.

**IMPORTANT!** Avoid creating bubbles during mixing and aliquoting.

- d. Select the program MagMAX\_Saliva\_2mL\_DUO on the instrument.
- e. Start the run, then load the prepared plates into position when prompted by the instrument.
- f. At the end of the run, immediately remove the plate from the instrument, then cover the plate or transfer the eluate to the final tube or plate of choice for final storage.

The purified DNA is ready for immediate use. Alternatively, store the plate at –20°C for long-term storage.

### Quantitation

To most accurately quantitate gDNA samples that are isolated from saliva, it is recommended to quantitate using either the Qubit<sup>™</sup> dsDNA BR (Broad Range) Assay Kit (Cat. No. Q32850) or Qubit<sup>™</sup> dsDNA HS (High Sensitivity) Assay Kit (Cat. No. Q32851). Another acceptable method is quantitation using qPCR and the Applied Biosystems<sup>™</sup> TaqMan<sup>®</sup> RNase P Detection Reagents Kit (Cat. No. 4316831).

#### Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



Manufacturer: Life Technologies Corporation | 2130 Woodward Street | Austin, TX 78744

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Revision history: Pub. No. MAN0017722

Revision	Date	Description
A.0	10 May 2018	New document

Important Licensing Information: This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.