

Maxwell 16 FFPE Tissue LEV DNA Purification Kit

Improve turnaround time for DNA extraction from FFPE samples

Achieve immediate time savings and enhance workflow with automation

The new Maxwell 16 FFPE Tissue Low Elution Volume (LEV) DNA Purification Kit enables automated extraction of genomic DNA from FFPE samples on the Maxwell 16 instrument. The Maxwell 16 FFPE Tissue LEV DNA Purification Kit allows you to extract DNA from up to 16 samples in approximately 30 minutes following an overnight proteinase-K digestion.

Automating nucleic acid extractions decreases hands-on bench time and reduces pipetting errors and repetitive motion injuries. The Maxwell 16 System speeds front-end processing, enabling you to perform additional molecular analyses without adding personnel. Maxwell 16 System is three times faster compared to manual methods and provides consistent high yields.

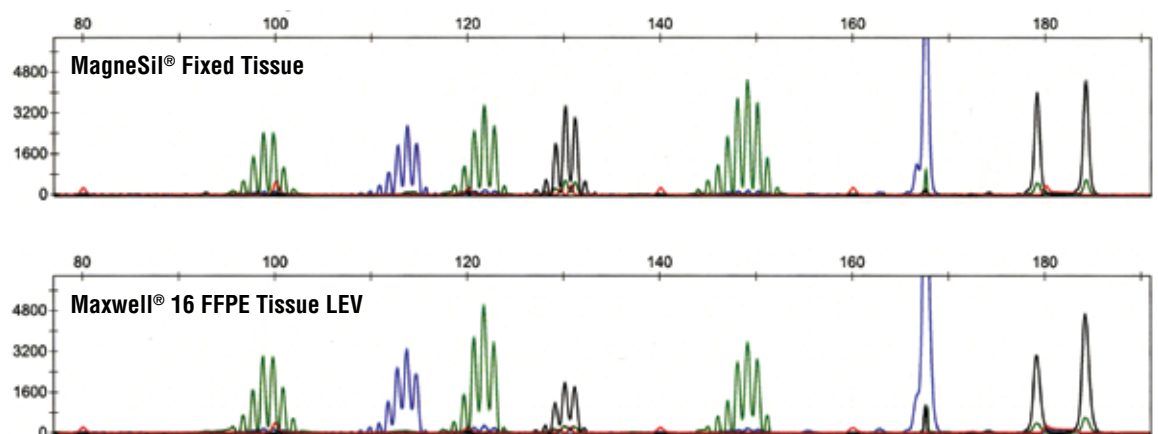


Figure 1. Gain confidence in your downstream analysis. DNA was purified from human FFPE tissue sections (5 micron) using MagneSil[®] Fixed Tissue System (A) or Maxwell[®] 16 FFPE Tissue LEV DNA Purification Kit (B) then analyzed using the MSI Analysis System, Version 1.2 (Cat. #MD1641) on an ABI 3100 capillary electrophoresis instrument.

Streamline purification for consistent results

The Maxwell 16 FFPE Tissue LEV DNA Purification Kit produces highly pure genomic DNA that is virtually free of PCR inhibitors without the use of harsh chemicals. Whether you are performing genetic screening assays such as Microsatellite Instability (MSI) Analysis, or correlating traditional morphology assessments with new genomics/genotyping techniques, the Maxwell 16 FFPE Tissue LEV DNA Purification Kit is a fast, safe and effective method for extracting biopsy specimens.

No cross-contamination

The Maxwell 16 FFPE Tissue LEV DNA Purification kit used in combination with the Maxwell 16 Instrument gives you the confidence that you can get consistent, high-quality results quickly. The unique design of the Maxwell 16 Instrument eliminates cross-contamination and common problems associated with other automated purification systems.

Maxwell[®] 16 FFPE Tissue LEV DNA Purification Kit

	Sample Avg. (pg/μl)	IPC (Ct)
Tissue A	49 ± 3	20
Tissue B	13 ± 8	20
Tissue C	3 ± 1	20

Get genomic DNA virtually free of PCR inhibitors. DNA was purified from FFPE sections (5 micron) from 3 different human tissues using the Maxwell[®] 16 FFPE Tissue LEV DNA Purification Kit then quantified using the real-time PCR Plexor[®] HY System (Cat.# DC1000). An internal PCR control (IPC) was used to test for the presence of PCR inhibitors. Data indicate none of the samples tested affected amplification of the IPC.

Visit www.maxwell16.com for more information.

Contact your Promega representative for product availability in your area.



7865CA

Maxwell® FFPE Tissue LEV DNA Purification Kit Specifications

Proteinase-K Digestion	Overnight
Processing Time	28 minutes
Extractions per Kit	48
Storage Requirements	22–25°C
Number of Samples per Run	Up to 16
LEV Configuration	20–100µl elution volume
General Purpose Laboratory Equipment	US only

Ordering Information

For specific countries contact your Promega representative

Product	Size	Cat#
Maxwell® 16 FFPE Tissue LEV DNA Purification Kit*	48 preps	AS1130
Maxwell® 16 Cell LEV DNA Purification Kit*	48 preps	AS1140
Maxwell® 16 Blood DNA Purification Kit*	48 preps	AS1010
Maxwell® 16 Cell DNA Purification Kit*	48 preps	AS1020
Maxwell® 16 Tissue DNA Purification Kit*	48 preps	AS1030
Maxwell® 16 LEV Cell Total RNA Purification Kit*	48 preps	AS1225
Maxwell® 16 LEV Tissue Total RNA Purification Kit*	48 preps	AS1220
Maxwell® 16 Total RNA Purification Kit*	48 preps	AS1050
Maxwell® 16 Instrument*	each	AS2000
Maxwell® 16 SEV Hardware Kit	each	AS1200
Maxwell® 16 LEV Hardware Kit	each	AS1250
Microsatellite Instability (MSI) Assay, Version 1.2	50 tests	MD1641

* For Laboratory Use



Promega Corporation • 2800 Woods Hollow Road • Madison, WI 53711-5399 USA • Telephone 608-274-4330 • Fax 608-277-2601
www.promega.com

©2008 Promega Corporation. All Rights Reserved.
 Prices and specifications subject to change without prior notice.

Printed in USA 11/08
 17056-FI-IN
 Part #1337

