

Maxwell 16 Cell LEV DNA Purification Kit

Streamline efficiency for your gDNA extraction from low cell volume samples

Consistent high yield from difficult samples

The Maxwell 16 Cell Low Elution Volume (LEV) DNA Purification Kit extracts genomic DNA from sample types inherently low in cell number including:

- Cerebral Spinal Fluid (CSF)
- Amniotic Fluid
- Cell Culture Supernatants

The system provides highly pure genomic DNA for downstream analysis. Additionally, the Maxwell 16 LEV format delivers a final DNA concentration that minimizes the need for additional processing of the extract prior to amplification. Consistent high yields and purity give you results you can trust. The Maxwell 16 System provides the easiest sample purification assuring confidence in your results and reducing the likelihood of time-consuming rework.

Immediately increase efficiency in your lab

The Maxwell 16 Cell LEV DNA Purification Kit automatically extracts DNA from low cell volume samples on the Maxwell 16 instrument. The system produces highly pure gDNA in about 30 minutes, instantly improving time to results.

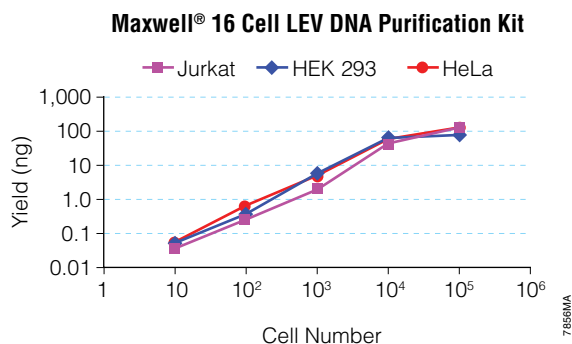
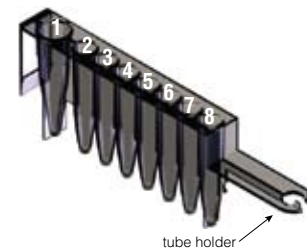


Figure 1. The Maxwell[®] 16 Cell LEV DNA Kit delivers consistent yields for amplification. DNA isolated from 10, 100, 1000, 10,000, and 100,000 cells provides yields required for downstream testing from as few as 10 cells. Purified DNA quantified using the Plexor[®] HY System (Cat. #DC1000).

Unique design eliminates cross-contamination

The Maxwell 16 System design eliminates cross-contamination and common problems associated with other automated purification systems, such as clogged tips and partial reagent transfers. Samples are placed directly into the cartridges and genomic DNA ready for amplification is obtained in approximately 30 minutes. Maxwell 16 uses a paramagnetic-particle handling system that does not require further liquid transference so there are no aerosols, drips or splashing.



Well Contents	User Adds:
1. Lysis Buffer	Sample
2. Paramagnetic Silica Particles	
3. Lysis Buffer	
4. Wash Buffer	
5. Wash Buffer	
6. Wash Buffer	
7. Empty	
8. Empty	Plunger

Figure 2. Maxwell[®] 16 LEV Cartridge

Visit www.maxwell16.com for more information.

Contact your Promega representative for product availability in your area.



Maxwell® Cell LEV DNA Purification Kit Specifications

Processing Time	30 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 Cell LEV DNA Purification Kit*	48 preps	AS1140
Maxwell® 16 FFPE Tissue LEV DNA Purification Kit*	48 preps	AS1130
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

*For Laboratory Use

