

Isolation of: **genomic DNA**

Sample Type: formalin-fixed paraffin-embedded (FFPE) tissue

Kit: Maxwell® 16 FFPE Plus LEV DNA Purification Kit **AS1135**

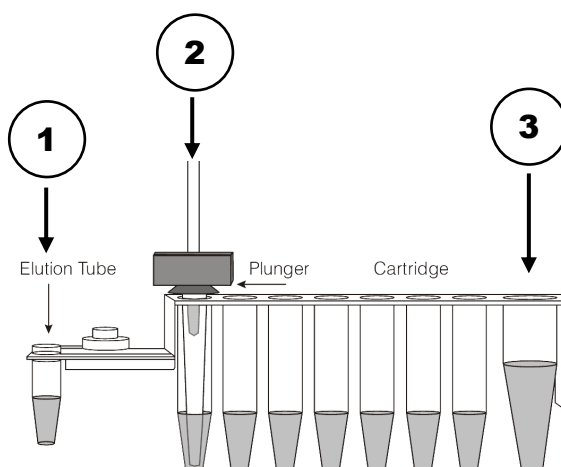
Optimized for
high yield. For
high purity use
AS1130.

Sample preparation:

- 1.) Dissolve supplied lyophilized Proteinase K by adding **500 µl** of nuclease-free water to each tube (final concentration 20 mg/ml). You can dispense the Proteinase K solution into smaller aliquots and store at –20°C for up to 1 year.
- 2.) Transfer your material (one to ten 5 µm sections from the FFPE sample) into a suitable microtube and overlay it with **180 µl** of the enclosed incubation buffer and **20 µl** of the Proteinase K solution from 1.). The sections should be completely covered by the solution!
- 3.) Incubate the sample at least 1 h at 70°C (higher yields are received by incubation for 8 h or overnight)
- 4.) Add **400 µl** of the included lysis buffer and vortex sample briefly.

Extraction:

- 1.) Place the cartridge to be used into the Maxwell LEV cartridge rack and remove the seal.
- 2.) Place one of the supplied elution tubes into the sample rack and add **50 µl** of the supplied elution buffer (①).
- 3.) Place the plunger in the indicated position of the cartridge (②).
- 4.) Transfer the complete sample into well 1 of the Maxwell cartridge (see image below ③).
- 5.) Select **LEV** configuration of the Maxwell and select method: **DNA → FFPE/Cells**. Start run.
- 6.) After the extraction your sample is ready-to-use for your downstream applications.



Further information can be found in the technical manual available online at
www.promega.com/resources/protocols