

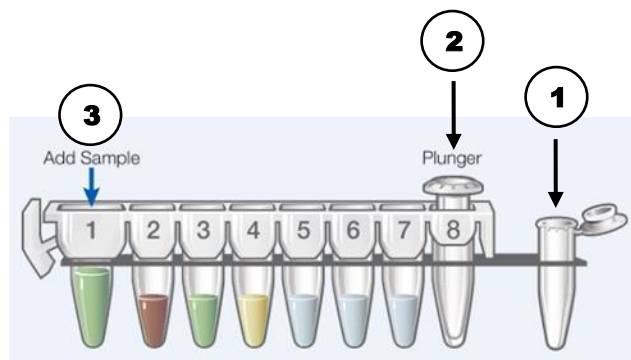
Isolation of: viral DNA
for detection of: HPV
Sample Type: swab (dry or ThinPreps)
Kit: Maxwell RSC Blood DNA Kit **AS1400**

Sample preparation:

- 1.) Add **400µl** water or NaCl to dry swab and incubate for 10 min. Carefully squeeze out liquid from swab in a 1.5ml tube.
If a swab with transport media is used, directly pipet **300 µl** of media to a 1.5ml tube.
- 2.) Add **300µl** lysis buffer and **30µl** Proteinase K.
- 3.) Vortex and incubate 20 min at 56°C.

Extraction:

- 1.) Place the cartridge to be used into sample rack and remove the seal.
- 2.) Place one of the supplied elution tubes into the sample rack and (see **①**).
- 3.) Place the plunger in the indicated position of the cartridge (see **②**).
- 4.) Add **60µl** of the supplied Elution Buffer into elution tube (see **①**).
- 5.) Transfer complete sample to the first well of the cartridge (see **③**).
- 6.) Select the appropriate program for the **Blood DNA Kit** and start run.
- 7.) After the extraction your sample is ready-to-use for your downstream applications.



Please note: This protocol is an instruction advice/recommendation but cannot replace in-house evaluation. Further information can be found in the technical manual available online at www.promega.com/resources/protocols