

Automated purification of *Aspergillus* DNA from blood samples

Aspergillus DNA was purified from blood samples using Maxwell® RSC Instrument

Kit:	Maxwell® RSC Blood DNA kit (Cat.# AS1400)
Analyses:	qPCR
Sample Type(s):	human whole blood
Input:	300µl
Materials Required:	<ul style="list-style-type: none">▪ Maxwell® RSC Instrument (Cat.# AS4500)▪ Maxwell® RSC Blood DNA Kit (Cat.# AS1400)▪ Lyticase▪ Thermoblock

This protocol was developed by Promega Applications Scientists and is intended for research use only.

Users are responsible for determining suitability of the protocol for their application.

For further information, see Technical Manual TM419, available at:

www.promega.com/protocols

or contact Technical Services at:
techserv@promega.com

Protocol:

1. Add 300µl of blood and 12.5µl of lyticase (4 U/µl) in a 1.5ml tube.
2. Incubate at 30°C and 1000rpm for 60 minutes.
3. Add 30µl of Proteinase K Solution to incubation tube.
4. Add 300µl of Lysis Buffer to incubation tube.
5. Vortex for 10 seconds.
6. Incubate at 56°C, 1000rpm for 20 minutes. Transfer each blood lysate sample from the incubation tube to well #1 of each cartridge. (Well #1 is the well closest to the printed side and furthest from the elution tube.)
7. Place a plunger in well #8 of each cartridge. Well #8 is the well closest to the elution tube.
8. Place an empty elution tube into the elution tube position for each cartridge. Add 50µl of Elution Buffer to the bottom of each elution tube.
9. Run the protocol RSC Blood DNA method on the Maxwell® RSC Instrument.

Results:

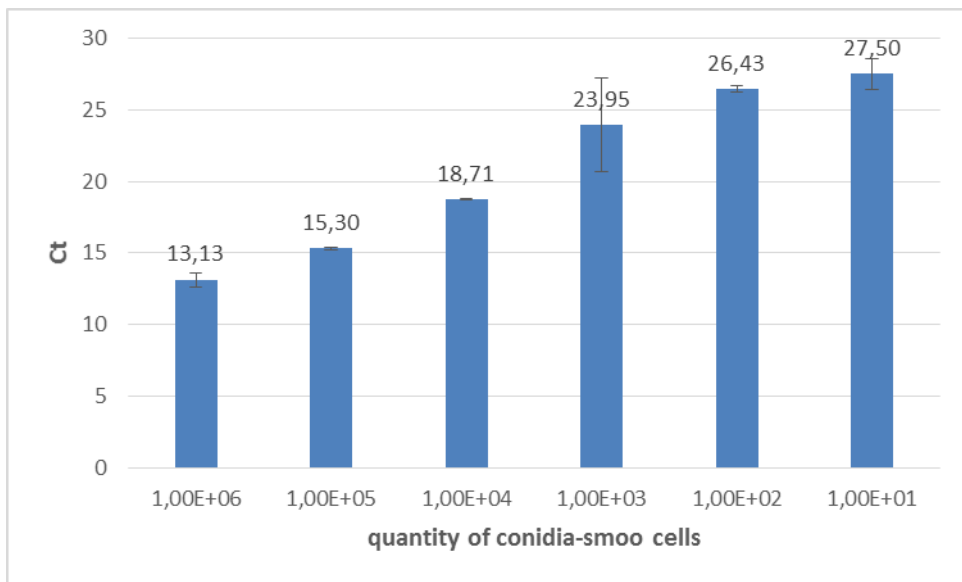


Figure 1. *Aspergillus* DNA detection in blood samples. Amplification of DNA extracted from 10^6 down to 10 conidia-smoo cells of *Aspergillus niger* spiked in blood samples using GoTaq® qPCR Master Mix (Cat.# A6001) and 28Sr specific *Aspergillus* primers¹. N=3. No amplification was observed in the non-template control.

Reference:

1. E. C. M. Williamson et al.(2000) PCR Diagnosis of Invasive Aspergillosis in BMT Recipients British Journal of Haematology, , 108, 132-139.