

KingFisher® Flex process with 96 deep well magnet head and NucleoMag™ 96 Blood Kit (Macherey-Nagel)

Sample process

1. Fill the plates according to the **Table 1**. More information about the reagents in the NucleoMag™ 96 Blood Kit instructions, Macherey-Nagel, catalog no. 744 500

Plate number	Plate type	Plate name	Content	Sample/reagent volume
1	Microtiter deep well 96 plate	Lysis	Lysis Buffer MB1 Blood	125 µl 100 µl
Dispense step, add...				
1			Binding Buffer MB2 NucleoMag B-Beads	360 µl 14 µl
2	Microtiter deep well 96 plate	Wash_MB3	Wash Buffer MB3	600 µl
3	Microtiter deep well 96 plate	Wash_MB4	Wash Buffer MB4	600 µl
4	Microtiter deep well 96 plate	Wash_MB5	Wash Buffer MB5	600 µl
6	KingFisher 96 plate	Elution	Elution Buffer MB6	100 µl

Table 1. Filling the plates

2. Combine the 96 deep well tip comb and the KingFisher 96 plate. For more detailed instructions, see KingFisher Flex User manual
3. Start the “**NucleoMag BloodDNA Flex96**” protocol using **arrow keys** and **START** button. You can also run the protocol using a computer, for more details see BindIt software user manual
4. Load the plates according to the protocol request and press **START** after every plate to confirm the action

Note! Confirm that the plates are placed in correct orientation: A1 well to be pointed to upper right corner of the plate holder in turntable. A1 row of the plate is then always located in the inner circle of the turntable
5. The purification protocol will start when the last plate is loaded and **START** button is pressed.
6. **Dispense step:** After lysis add 14 µl of resuspended NucleoMag B-Beads and 360 µl of Binding Buffer MB2 to plate 1. Beads and MB2 may be premixed before addition to plate 1
7. After the purification process is completed the plates are removed according to instructions shown in the instrument screen. Press **START** after each plate removal to confirm the action
8. When the last plate is removed text End of run will appear. Press **STOP** to complete the run.

For more information: www.thermoscientific.com/kingfisher and www.mn-net.com