

KingFisher® Flex process with 24 magnet head and InviMag® Blood DNA Kit (Stratec Molecular)

Sample process

1. Fill the plates according to the **Table 1**. More information about the reagents in the InviMag® Blood DNA Kit Manual, Stratec Molecular, catalog no. 74313001.

Plate number	Plate type	Plate name	Content	Sample/reagent volume
1	KingFisher Flex 24 deep well plate	Lysis / Binding	Lysis Buffer A Blood Proteinase K	1000 µl 1000 µl 100 µl
Dispense step, add...				
1			Binding buffer B6 MAP Solution A	2000 µl 80 µl
2	KingFisher Flex 24 deep well plate	Wash 1	Wash Buffer I	2500 µl
3	KingFisher Flex 24 deep well plate	Wash 2	Wash Buffer I	2500 µl
4	KingFisher Flex 24 deep well plate	Wash 3	Wash Buffer II	4000 µl
5	KingFisher Flex 24 deep well plate	Wash 4	Wash Buffer II	4000 µl
6	KingFisher Flex 24 deep well plate	Elution	Elution Buffer D	500 µl

Table 1. Filling the plates

2. Start the “**InviMag_BloodDNA_Flex24**” protocol using **arrow keys** and **START** button. You can also run the protocol using a computer, for more details see BindIt software user manual.
3. Load the plates according to the protocol request and press **START** after every plate to confirm the action. Use KingFisher Flex 24 deep well tip comb and plate for tip loading.

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.

4. The purification protocol will start when the last plate is loaded and **START** button is pressed.
5. **Dispense step:** After lysis add 80 µl of resuspended MAP solution A and 2000 µl of Binding Buffer B6 to plate 1. Beads and Binding buffer may be premixed before addition to plate 1.
6. 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.
7. 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.invitek.de