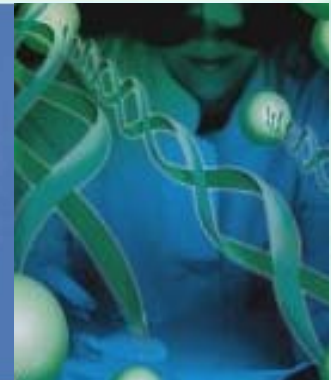




Gene expression studies with MagRo™ 96-M

Automated solutions for
RNA & mRNA purification



Bio-Nobile's PickPen magnetic particle technology:

- saves time effort as elution step is eliminated
- 100% transfer of all purified mRNA into RT-PCR

The MagRo 96-M workstation quickly and efficiently processes all 96 samples in a standard or deep well plates. Using the proven PickPen® method the instrument's fast protocols allow as many as 2000 purifications a day. Open design allows for easy integration into laboratory robotics.



[Learn more about PickPen technology](#)

[Learn more about MagRo 96-M](#)



After purification reporter gene assays can be easily performed with the BERTHOLD TECHNOLOGIES multimode and dedicated microplate readers.

[Learn more about microplate readers](#)

[Get in touch with Berthold Technologies](#)



Visit us at the Biotechnica in Hannover, Germany:
BERTHOLD TECHNOLOGIES Hall 9, Booth E02
BIONOBILE Hall 9, Booth E010/1