

## Can MicroPlex technology detect mRNA?

It is possible to detect mRNA by using RT-PCR amplified material. Direct detection of mRNA may be possible with an appropriate assay design.

Luminex Technical Applications Group has detected PCR-amplified cDNA at 15 and 30 cycles of amplification when using 20 ng and 4 ng of cDNA template, respectively. Median Fluorescent Intensity (MFI) greater than two times the MFI of the background (no sample) wells was considered positive.

