

PRODUCT NAME		
Human ChIP-seq grade Myoglobin Exon 2 primer pair		
Cat. No: <b>pp-1006-050</b>	Format: 50 $\mu$ l	Concentration: 10 $\mu$ M
Cat. No: <b>pp-1006-500</b>	Format: 500 $\mu$ l	Concentration: 10 $\mu$ M

**Product Description:** This primer pair specifically amplifies a genomic region from exon 2 of the human myoglobin gene. These Primers are thoroughly tested and optimized for routine SYBR® Green Real-Time qPCR assay following ChIP and for ChIP-sequencing library validation (e.g. before and after ChIP-seq library preparation).

**Amplicon length:** 76 base pairs.

**Amplified region:** chr22:34,336,993-34,337,068

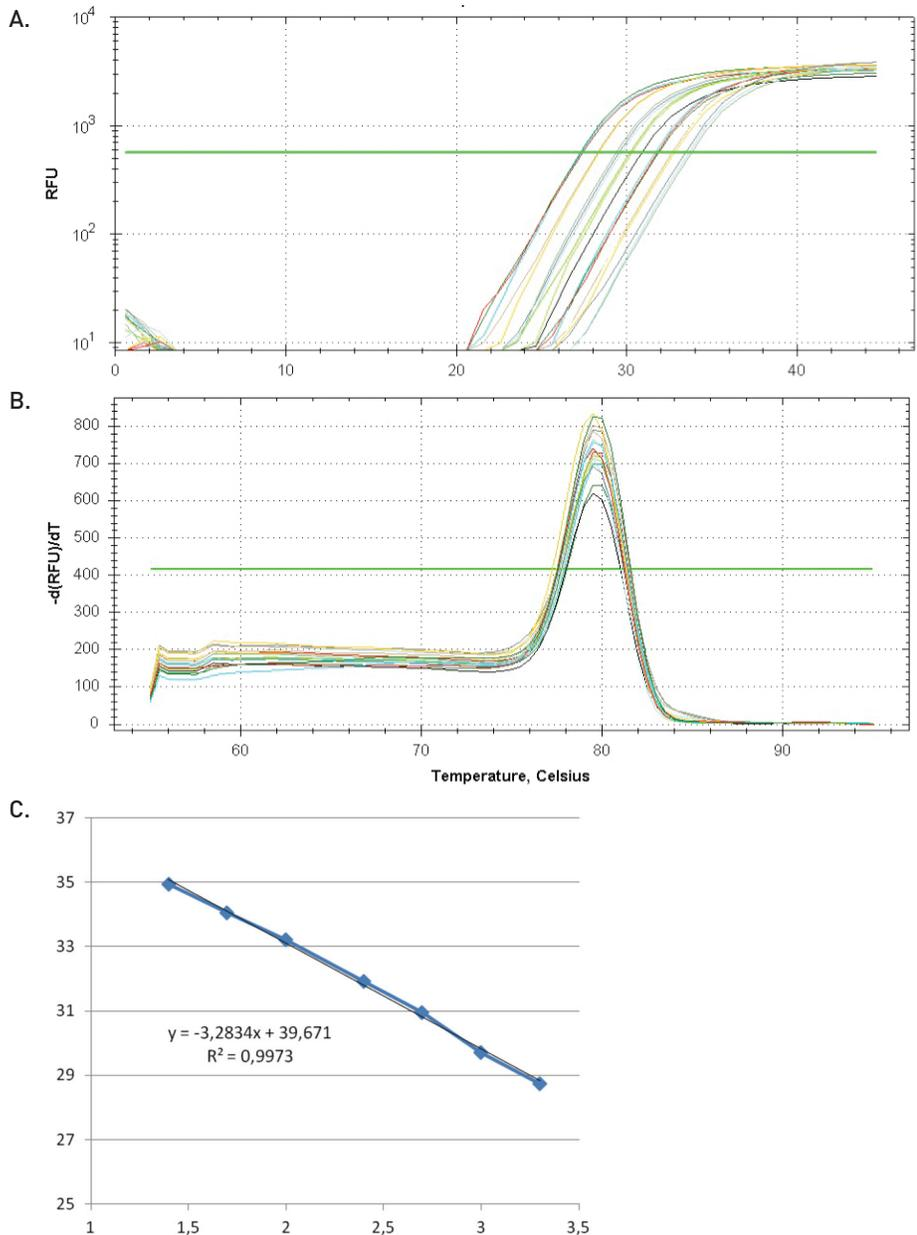
**Specificity:** Human

**Format:** 10  $\mu$ M solution in MilliQ water (5  $\mu$ M of each primer)

**Storage:** For long storage, store at -20°C. Avoid multiple freeze-thaw cycles.

**Precautions:** This product is for research use only. Not for use in diagnostic or therapeutic procedures.

**Last data sheet update:** April 18, 2013



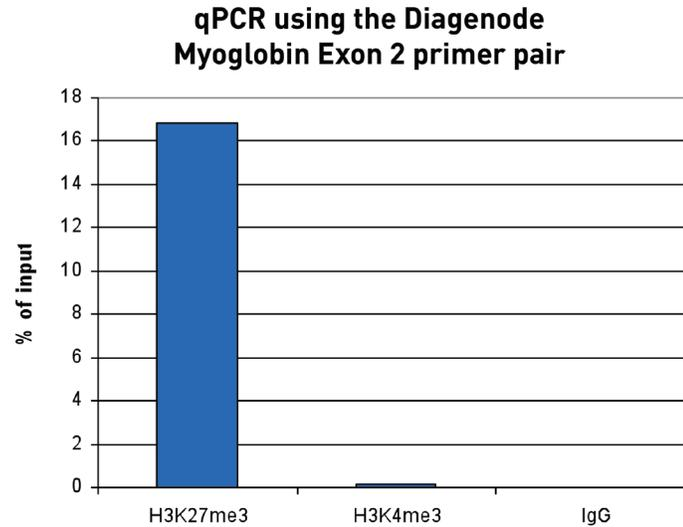
**Figure 1**

Sheared DNA from HeLa cells was analysed in triplicate by SYBR® Green real-time PCR using the Diagenode MB ex2 Primer Pair [cat. No. pp-1006-050, pp-1006-500]. A dilution series ranging from 2 ng to 25 µg of DNA template was amplified with 1 µl of the provided primers in 25 µl total reaction volume on an iQ5 thermocycler (Biorad). qPCR conditions were as follows: incubation at 95°C for 10 minutes, followed by 45 cycles of 30 seconds at 95°C, 30 seconds at 60°C, and 30 seconds at 72°C and a final extension for 10 minutes at 72°C.

Figure 1A: amplification curves [logarithmic view]. The green line represents the position of the threshold.

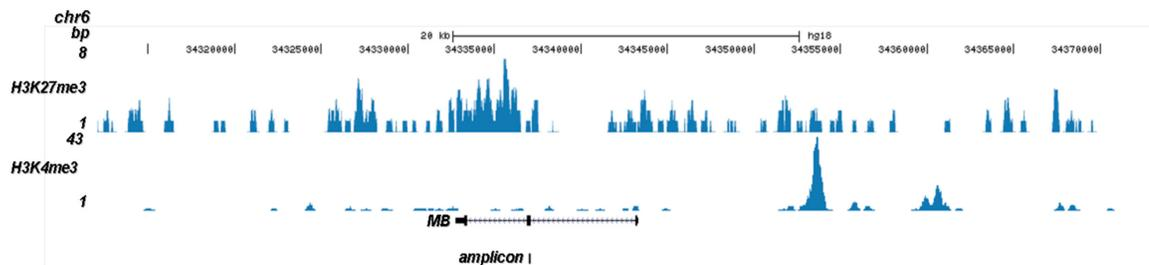
Figure 1B: melting curve analysis of the different amplification products.

Figure 1C: plot of the Ct value [mean of 3 replicates] against the log of the initial DNA amount. The reaction efficiency, as calculated from the slope of this curve, is 101.6%.



**Figure 2**

ChIP was performed on HeLa cells using antibodies against H3K27me3 (cat. No. pAb-069-050), located at inactive genomic regions, and H3K4me3 (cat. No. pAb-003-050), known to be located at active promoters. Rabbit IgG was used as a negative IP control. The ChIP'd samples were analysed by qPCR using the MB ex2 Primer Pair (cat. No. pp-1006-050, pp-1006-500). Figure 2 shows the recovery, expressed as a % of input (the relative amount of immunoprecipitated DNA compared to input DNA after qPCR analysis).



**Figure 3**

ChIP was performed as described above and the ChIP'd DNA was subsequently analysed by high throughput sequencing on an Illumina GAIIx. Figure 3 shows the H3K27me3 and H3K4me3 profiles in a region of chromosome 6 containing the MB gene. The position of amplicon obtained with the Diagenode MB ex2 Primer Pair (cat. No. pp-1006-050, pp-1006-500). is indicated.