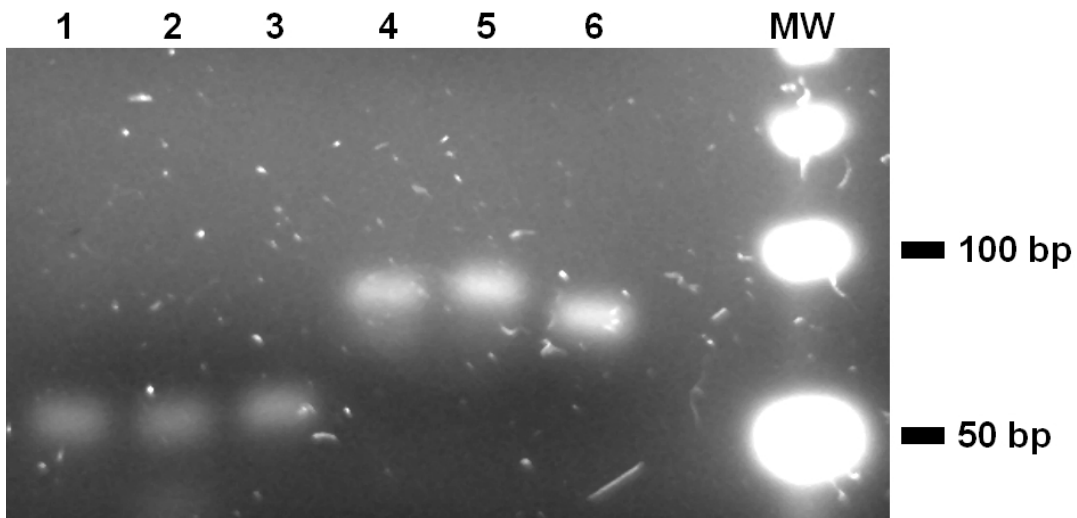


**Figure 1:** DNA from undifferentiated human teratocarcinoma NCCIT cells was analyzed in duplicate by real-time PCR starting from 5  $\mu$ l of DNA template (0.01 $\mu$ g/ml) using the Diagenode primers to amplify a region located at -1.0kb of the human GAPDH promoter (cat#: pp-1003-050, -500). One  $\mu$ l of provided primer pairs is used by PCR of 25  $\mu$ l final volume. A Real-Time PCR Detection System and iQ SYBR Green have been used. qPCR conditions used are as follows: 95°C for 3 minutes, 40 cycles of: [95°C for 15 seconds, 60°C for 45 seconds] and 1 cycle of 95°C for 1 minute. Duplicates are shown in blue. Threshold position is in orange.



**Figure 2:** qPCR products were analysed by electrophoresis (2% agarose gel) stained with SYBR Safe and illuminated with UV light. The right lane shows molecular weight markers (MW) that decrease in size by 50 bp. Different qPCR products using different primer pairs which are available at Diagenode were tested: 1: primers for human GAPDH promoter (cat#: pp-1001-050 and cat#: pp-1001-500), 2: primers for human GAPDH promoter -0.5kb (cat#: pp-1002-050 and cat#: pp-1002-500), **3: primers for human GAPDH promoter -1.0kb (cat#: pp-1003-050 and cat#: pp-1003-500)**, 4: primers for human c-fos promoter (cat#: pp-1004-050 and cat#: pp-1004-500), 5: primers for human beta-actin promoter (cat#: pp-1005-050 and cat#: pp-1005-500) and 6: primers for human myoglobin exon 2 (cat#: pp-1006-050 and cat#: pp-1006-500).