A C T I V E 🚺 M O T I F®

ChIP-IT[®] Control Kit – Mouse

Catalog No: 53011 Format: 5 rxns

Quality Control: ChIP-IT[®] Control Kit – Mouse is quality control tested in combination with Active Motif's ChIP-IT[®] Express Kit (Catalog No. 53008).

Mouse macrophage 4/4 cells were grown, fixed and used to prepare chromatin as described in the ChIP-IT Express manual. ChIP reactions were then performed using 2 μ g RNA pol II antibody plus 2 μ g bridging antibody or 2 μ g negative control IgG. The immunoprecipitated DNA and the control Input DNA were then used in endpoint PCR using the EF1-alpha control primers (Figure 1). The reactions were cycled for 36 repetitions. The positive control primers generate a 233 bp product which should be enriched in the RNA pol II and Input samples. Signal in the Negative IgG samples represents non-specific background.

Kit Components:

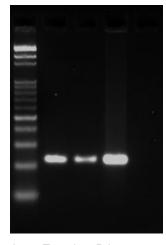
50 μl RNA pol II mouse monoclonal antibody (0.2 μg/μl) (also sold as Cat. No. 39097)
50 μl Bridging antibody (1 μg/μl) (also sold as Cat. No. 53017)
50 μl Negative control mouse IgG (0.2 μg/μl)
400 μl EF1-alpha primer mix (2.5 μM)
1.5 ml 10X PCR buffer
1.5 ml 10X PCR loading dye

Endpoint PCR Analysis We recommend the following PCR conditions: 9.8 μl DEPC H₂O 2.5 μl 10X PCR Buffer 2.5 μl 10X PCR Loading Dye 1.0 μl dNTPs (5 mM mix) 0.2 μl Taq polymerase 4.0 μl EF1-alpha primer mix 5.0 μl ChIP DNA 25 μl Total Volume

Reactions were cycled 36 times with the following steps per cycle: 94°C denaturing for 20 seconds 59°C annealing for 30 seconds 72°C extension for 30 seconds

Storage and Guarantee: The ChIP-IT Control Kit – Mouse components are shipped on dry ice. The negative control IgG antibody should be stored at 4°C, all other components can be stored at -20°C.

This product is guaranteed for 6 months from date of receipt under the correct storage conditions. Aliquot antibodies to avoid exposing to multiple freeze-thaw cycles.



Lane Template Primers 1 DNA Ladder --2 RNA pol II EF1-alpha 3 Negative IgG EF1-alpha 4 Input DNA EF1-alpha 5 H₂O control EF1-alpha