## ChIP-IT® Control qPCR Kit – Rat



Catalog No: 53028

Format: 5 rxns

**Quality Control:** ChIP-IT<sup>®</sup> Control qPCR Kit – Rat is quality control tested in combination with Active Motif's ChIP-IT<sup>®</sup> Express Kit (Catalog No. 53008).

Chromatin from a rat sample is fixed and sheared as described in the ChIP-IT Express manual. 5  $\mu$ g of the prepared chromatin is used for each ChIP reaction. Positive control ChIP reactions are performed in triplicate using 2  $\mu$ g RNA pol II antibody plus 2  $\mu$ g Bridging antibody, while negative control ChIP reactions are performed in triplicate using 2  $\mu$ g negative control antibody. The immunoprecipitated DNA and the Input DNA are used in qPCR with both Actb and Negative-1 qPCR primer sets.

To pass QC, the RNA pol II ChIP reactions should produce at least 5-fold enrichment with the Actb primer set as compared to the Negative-1 primer set. The negative control ChIP reaction should not show enrichment greater than half the RNA pol II enrichment with the Actb primer set (Figure 1).

## Contents:

RNA pol II mouse monoclonal antibody (0.2  $\mu$ g/ $\mu$ I) (also sold as Catalog No. 39097) Bridging antibody (1  $\mu$ g/ $\mu$ I) (also sold as Catalog No. 53017) Negative control mouse IgG (0.2  $\mu$ g/ $\mu$ I) Actb qPCR primer set (2.5  $\mu$ M) (also sold as Catalog No. 71026) Negative-1 qPCR primer set (2.5  $\mu$ M) (also sold as Catalog No. 71024)

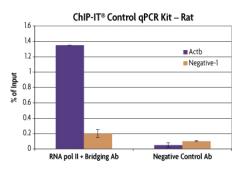
## qPCR Analysis

We recommend the following qPCR conditions as a starting point. Please optimize for your qPCR machine and master mix as needed.

SYBR green Master Mix 5 µl
PCR primer mix (2.5 µM) 0.7 µl
Sterile water 1.3 µl
ChIP DNA 3 µl
Total Volume 10 µl

95°C for 2 min 40 cycles as follows: 95°C for 15 sec 58°C for 20 sec 72°C for 20 sec

**Storage and Guarantee:** The ChIP-IT® Control qPCR Kit – Rat 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 the antibodies to avoid exposing to multiple freeze-thaw cycles.



**Figure 1:** qPCR analysis of 5 µg rat brain chromatin assayed using the ChIP-IT Control qPCR Kit – Rat and ChIP-IT Express Kit.