

RNA ChIP-IT[®] Control Kit – Human

Catalog No: 53025

Format: 5 rxns

Quality Control: The RNA ChIP-IT Control Kit – Human is quality-control tested to demonstrate that the positive and negative control antibodies are effective in performing RNA-chromatin immunoprecipitation.

HeLa cells were grown, fixed and used to prepare chromatin as described in the RNA ChIP-IT Kit manual. RNA-ChIP reactions were then performed using 10 μ l of Suz12 antibody and 10 μ l of control rabbit IgG (0.2 μ g/ μ l). The immunoprecipitated RNA and the control input RNA were then used in RT-PCR using lincSFPQ primer mix.

Kit Components:

50 μ l Suz12 rabbit polyclonal antibody

100 μ l rabbit IgG (0.2 μ g/ μ l)

400 μ l lincSFPQ primer mix

Storage and Guarantee: The RNA-ChIP Control Kit – Human components are shipped on dry ice. The lincSFPQ primer mix, the Suz12 antibody and the control rabbit IgG should 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.

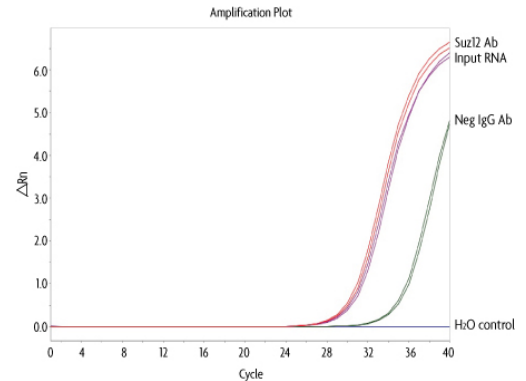


Figure 1: Amplification plot corresponding to qRT-PCR amplification of lincSFPQ locus from Suz12 antibody (Cat. No. 39357) and rabbit IgG RNA-ChIP samples.

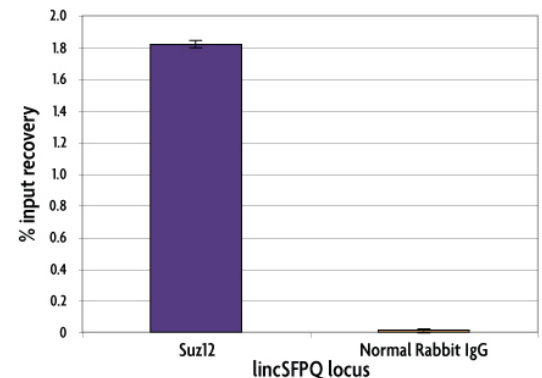


Figure 2: RNA-ChIP on HeLa chromatin using Suz12 antibody (Cat. No. 39357) and rabbit IgG. qRT-PCR is performed using primers amplifying lincSFPQ RNA locus. Results are shown in terms of percentage input recovery.