

## *Drosophila* Negative Control Primer Set 3

**Catalog No:** 71038

**Format:** 96 rxns

**Background:** The *Drosophila* Negative Control Primer Set 3 amplifies a 65 base pair fragment from 13 kb 3' of the start site of *D. melanogaster* split ends (Spen) transcript on chromosome 2L. It can be used as a negative control for spike-in experiments with H3K27me3 and EZH2.

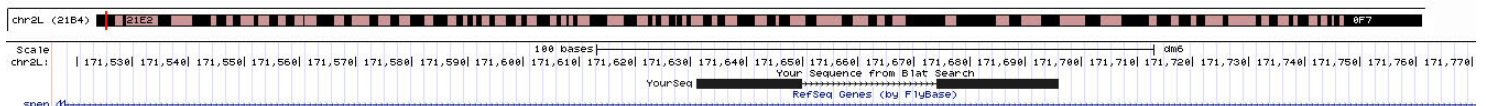
**Contents:** This control primer set contains both forward and reverse primers in 400 µl of nuclease-free distilled water. The final concentration for each primer is 2.5 µM.

**Application Notes:** Amplification should be carried out in a total volume of 20 µl, using the DNA template, 4 µl of the primer set, and 10 µl SYBR Green 2X qPCR Master mix with an annealing temperature of 58°C. For genomic DNA amplification, 12.5 ng of DNA was used as template.

**Quality Control:** This primer set was used to produce a single PCR product from genomic DNA using qPCR to generate an amplification curve with a Ct of fewer than 28 cycles. After amplification, melt curve analysis was performed to confirm the production of a single PCR amplicon.

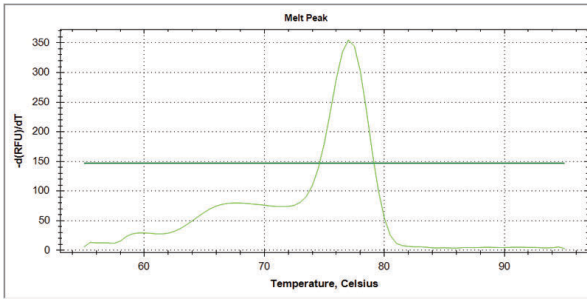
**Storage and Guarantee:** The primers are shipped at room temperature and should be stored at -20°C upon receipt to ensure stability. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



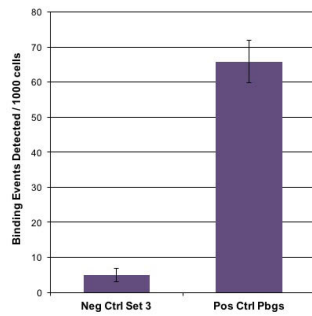
### Genomic Location

Image representing the relative location of the primer set amplicon within the genome, as generated by the UCSC Genome Browser.



### Melt Curve

PCR product melt curve for qPCR reaction using 12.5 ng of total DNA as template. The single peak corresponds to a single amplicon.



### ChIP qPCR Data

ChIP was performed on chromatin from *D. melanogaster* adults using an antibody to H3K27me3 and subjected to qPCR with the indicated primer set (X-axis). Data presented are normalized binding events per 1000 cells.