

## Zebrafish Negative Control Primer Set 1

**Catalog No:** 71034

**Format:** 96 rxns

**Background:** The Zebrafish Negative Control Primer Set 1 amplifies a 64 base pair fragment from a gene desert on zebrafish chromosome 3. It can be used as a control for almost all transcription factors, RNA pol II and most histone modifications.

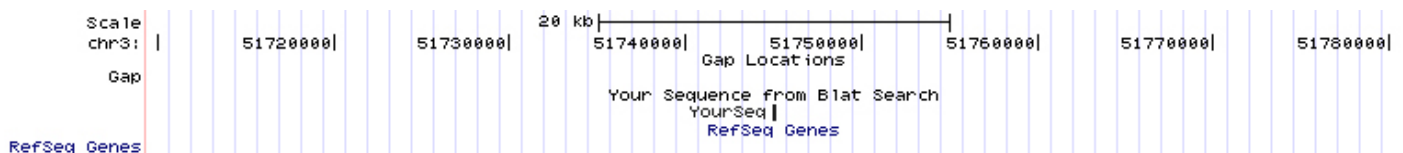
**Contents:** This control primer set contains both forward and reverse primers in 400 µl of nuclease-free distilled water. The final concentration for each primer is 2.5 µM.

**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.

