

## MQPROSHR-3 Datasheet

### Primer Information

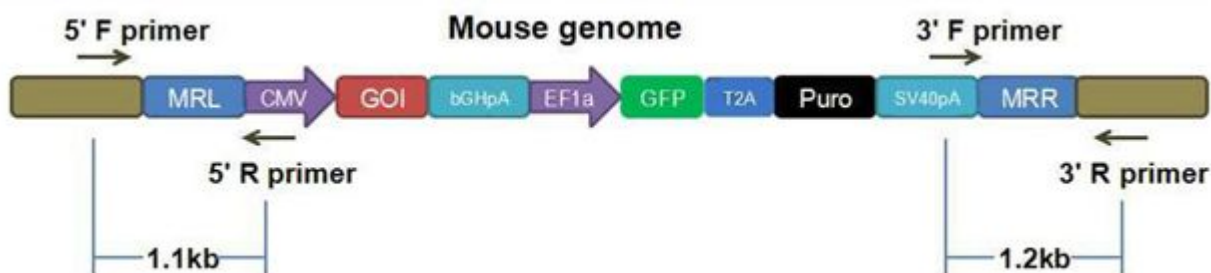
---

**Catalog No.:** MQPROSHR-3

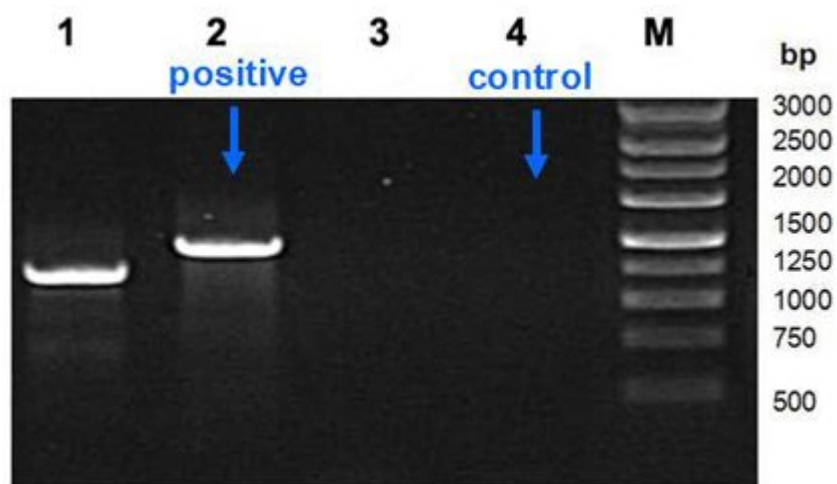
**Description:** 3' Junction-PCR Primer Set

**Forward Primer:** 5'-GGTGCCTGAAATCAACCTCTGGA-3'

**Reverse Primer:** 5'-TCAAGCCAGTCCAAGAGAAAGCA-3'



Note: PCR primers designed to amplify the HR junction were used to verify the successful integration.



Note: Positive (Lane 2) and negative control (Lane 4) PCR products were run on the 1% Agarose/EtBr gel.

---