

HQPAVSHR-5 Datasheet

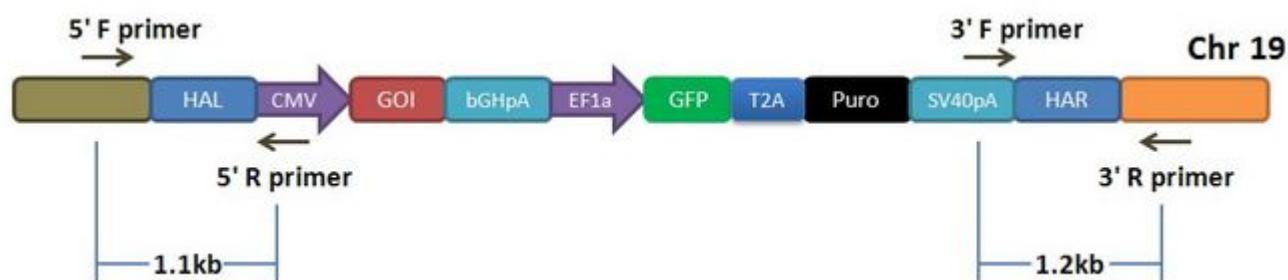
Primer Information

Catalog No.: HQPAVSHR-5

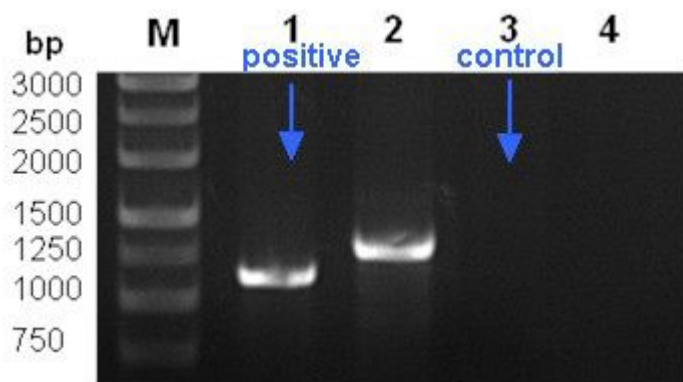
Description: 5' Junction-PCR Primer Set

Forward Primer: 5'-CCGGAAGTCTGCCCTCTAAC-3'

Reverse Primer: 5'-CCCGTGAGTCAAACCGCTAT-3'



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



Note: Positive (Lane 1) and negative control (Lane 3) PCR products were run on the 1% Agarose/EtBr gel.
