

SNAP-CHIP® OncoStat™ Full Panel Primer Set



EpiCypher®

Catalog No. 18-6201

Lot No.

Pack Size 100 reactions

Product Description:

The SNAP-CHIP® OncoStat™ Full Panel Primer Set contains a universal forward primer pre-mixed with reverse primers that uniquely amplify DNA-barcoded nucleosomes in the SNAP-CHIP OncoStat Panel (Catalog No. 19-2001), consisting of H3.3 wild-type (WT), H3.3K4M, H3.3K9M, H3.3K27M, H3.3G34R, H3.3G34V, H3.3G34W, and H3.3K36M. These primers can be used in SNAP-CHIP experiments where quantitative PCR (qPCR) provides an assessment of antibody performance and technical variability.

The primers are compatible with SYBR™ chemistry. For TaqMan® applications, the universal SNAP-CHIP Dual Label Hydrolysis Probe can be purchased separately (Catalog No. 18-6001). The volume is sufficient to perform 100 x 10 µL qPCR reactions (see SNAP-CHIP manual).

Formulation:

K-MetStat universal forward primer pre-mixed with individual reverse primers at 5 µM (20X) individual primer concentration in Molecular Grade H₂O.

Storage and Stability:

Stable for 2 years at -20°C from date of receipt.

Application Notes:

For use in SNAP-CHIP-qPCR. Compatible with the SNAP-CHIP OncoStat Panel (Catalog No. 19-2001) and the SNAP-CHIP Dual Label Hydrolysis Probe (Catalog No. 18-6001).

DNA Sequence:

SNAP-CHIP OncoStat Forward Primer 5'-CGTATCGCGCGCATAATA-3'

H3.3 WT Reverse Primer 5'-ATACGACGAGATAGTCGACG-3'

H3.3K4M Reverse Primer 5'-TCGCGCGTACGAAAC-3'

H3.3K9M Reverse Primer 5'-ATTCGCGCGTACGTATAC-3'

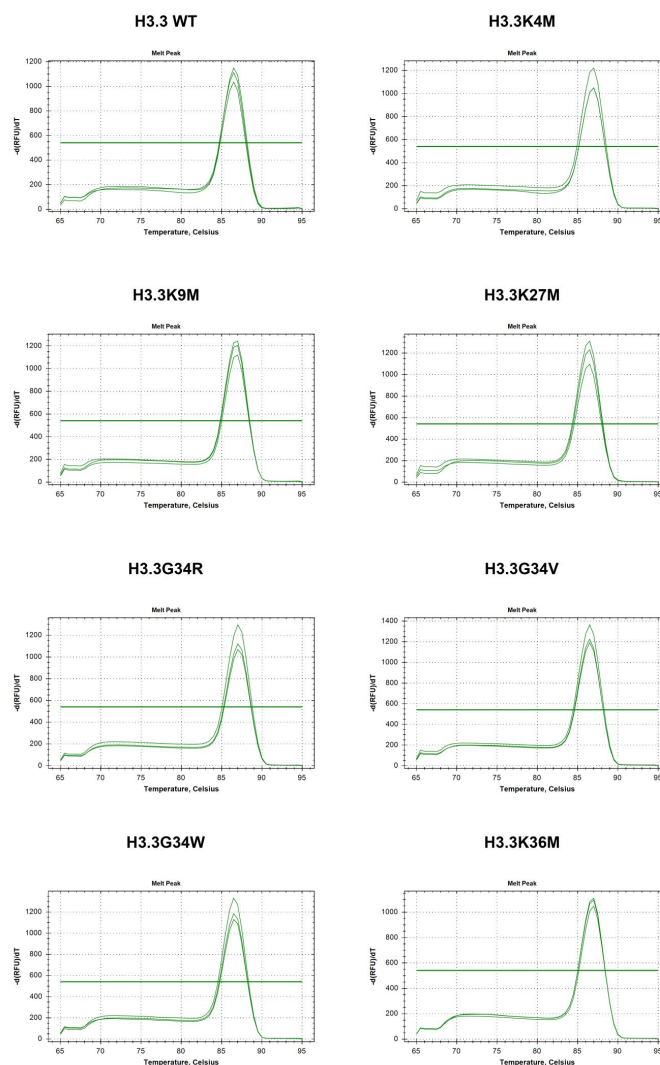
H3.3K27M Reverse Primer 5'-ACGCGATAACGTCGACTA-3'

H3.3G34R Reverse Primer 5'-TATCGCGTCGTACGATCG-3'

H3.3G34V Reverse Primer 5'-CGCGATTCGCGTAATACG-3'

H3.3G34W Reverse Primer 5'-GCGACGCGTAATCGA-3'

H3.3K36M Reverse Primer 5'-TAATCGACGCGTTACGC-3'



Melt curve: SYBR Green quantitative PCR using Input DNA from a chromatin immunoprecipitation experiment (ChIP-qPCR) where K-562 cell human chromatin was spiked-in with the SNAP-CHIP OncoStat Panel (EpiCypher Catalog No. 19-2001). PCR products of the indicated primer sets were subjected to a melting curve analysis, where the presence of a single peak indicates specific amplification of one product. Data shown is for triplicate qPCR reactions in a single ChIP experiment.

This product is for *in vitro* research use only and is not intended for use in humans or animals.