



Universal methylated DNA standard (mouse) Protocol

Catalog number: ENZ-45006

Primer Set

mMLH1 Primer F:

5' - GGTGTACGAAGTTATTTTTATTTTAGTC - 3'

mMLH1 Primer R:

5' - ACCCAACGATACCTAATAATAAAACC - 3'

PCR Setup:

The following setup is designed for a 25 μ l total reaction volume:

Component	Volume	Final Conc.
mMLH1 primer F*	Variable	0.2 to 0.8 μ M
mMLH1 primer R*	Variable	0.2 to 0.8 μ M
Bisulfite-converted DNA	2 μ l	up to 20 ng/ μ l
10 mM dNTP mix	0.5 μ l	0.2 mM each dNTP
Standard PCR buffer	Variable	1x
MgCl ₂ or MgSO ₄	Variable	1-4 mM, if needed
Hot-Start DNA Polymerase	1 to 2 units	
<i>Hot-start DNA polymerase)</i>	Variable 1 to 2 units	
Add water to 25 μ l		

* Alternatively, you may substitute primers of your choice.

** Remember to bisulfite-treat the DNA prior to performing PCR.

Recommended Thermocycler Conditions:

95 °C, 10 minutes

95 °C, 30 seconds

58 °C, 30 to 60 seconds

72 °C, 60 seconds

Repeat steps B through D an additional 29 to 39 times depending on the polymerase used.

72 °C, 7 minutes

4 °C

The PCR amplicon can now be used directly for sequencing analysis

Additional information including expected results

The expected PCR amplicon for the Universal Methylated Mouse DNA Standard is 304 bp, corresponding to nucleotide positions 430 to 778 of mouse MLH1 DNA including the regions (italicized) that hybridize to the primers (GenBank Accession #: AF400617).

Original sequence of mouse MLH1 DNA for bisulfite treatment and PCR amplification (sense strand 5' to 3'). The cytosines (underlined) in the CpG dinucleotide context (bold capital letters) are methylated enzymatically by M.Sss1 methyltransferase:

```

421 -----g gtgtaCGaag tcaccctcac cccagcCGCG
461 acccttcaag gccaagaagC GgcagaggC Gaggcctgcc
501 CGCGtCGctc tctcctCGg agtgagcaCG gCGgccaaag
541 acatgtcacc ctgcCGcaga CGctCGacca gggcCGCGCG
581 ttctCGtcc cctacaaacC GctCGtagaa ttCGtgctCG
621 gcctCGtagt ggCGcctcaC GtCGCGttcc CGagtagagg
661 CGaccaggCG gCGacacacc aggcacaggg cccCGtcacc
701 ctcCGcaggc tccaccacca ggtatCGctg ggt-----

```

Expected sequence of the above DNA following bisulfite treatment. Methylated cytosines in the CpG dinucleotide context remain unconverted following bisulfite treatment, whereas non-methylated cytosines, or cytosines not in the CpG context, are converted to uracil and detected as thymine after PCR.

```

421 -----g gtgtaCGaag ttatttttat ttagtCGCG
461 Atttttaag gtaagaagC GgtagaggtC Gaggttgtt
501 CGCGtCGttt ttttttCGg agtgagtaCG gCGgttaaag
541 atatgttatt ttgtCGtaga CGttCGatta gggCGCGCG
581 ttttCGttt ttataaatC GttCGtagaa ttCGtgttCG
621 gtttCGtagt ggCGttttaC GtCGCGtttt CGagtagagg
661 CGattaggCG gCGatatatt aggtataggg tttCGttatt
701 tttCGtaggt ttattatta ggtatCGttg ggt-----

```