

# Universal methylated DNA standard (mouse)

## Methylation standard (mouse) to help determine efficiency of a bisulfite DNA conversion

The Universal mouse methylated DNA standard with the included primers can be used in conjunction with Enzo's Blood and tissue methylation kit or the Express methylation kit. The DNA included in this set is enzymatically methylated at all cytosine positions to help assess the efficiency of the bisulfite-mediated conversion of DNA. DNA is isolated from a male mouse source. The primer set is designed to amplify a fragment of the mouse mismatch MLH1 mismatch repair gene following bisulfite treatment. The methylated cytosines comprising CG dinucleotides remain unconverted following bisulfite treatment, whereas non-methylated cytosines are converted into uracil and detected as thymine after PCR.

### Ordering Information

[Order Online »](#)

ENZ-45006-0001	20 tests
----------------	----------

### Manuals, SDS & CofA

[View Online »](#)

- Conveniently use with methylation kits
- Enzymatically modified standard includes short, easy to follow protocol
- Primers included

mMLH1 Primer I:

5' - GGTGTACGAAGTTATTTTATTTTAGTC - 3'

mMLH1 Primer II:

5' - ACCCAACGATACCTAATAATAAAACC - 3'

The following primer set is provided in the pack.  
Following bisulfite treatment of experimental methylated DNA, PCR can be conducted using the mMLH1 set primers with the universal methylated DNA as a control.

## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. With proper storage, good for one year upon receipt.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Contents</b>	Control primers: (mMLH1 (F) and mMLH1 (R)) 20µl each and Universal methylated DNA (mouse) 5µg/20µl
<b>Formulation</b>	Universal methylated DNA standard (mouse), 250ng/µL in TE buffer, pH 8.0. Control primers 20µM each primer in TE buffer, pH 8.0



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)