

## **BioArray™ 5-hmC Methylation Kit**

Catalog #: ENZ-45011

## **USE FOR RESEARCH PURPOSES ONLY**

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

## **LIMITED WARRANTY; DISCLAIMER OF WARRANTIES**

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NON-INFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

## **TRADEMARKS AND PATENTS**

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending. Enzo is a trademark of Enzo Life Sciences, Inc.

**FOR RESEARCH USE ONLY.  
NOT FOR USE IN DIAGNOSTIC PROCEDURES.**



Please read entire booklet before proceeding with the assay.



Carefully note the handling and storage conditions of each kit component.



Please contact Enzo Life Sciences Technical Support if necessary.

## TABLE OF CONTENTS

Introduction .....	2
Assay Principal .....	2
Materials Supplied .....	5
Storage .....	5
Precautions.....	5
Reagent Preparation .....	6
Methods & Procedures .....	6
Protocol .....	7
Appendix I:	
5-hmC Control DNA .....	10
Appendix II:	
Glucosylation of 5-hmC in DNA .....	11
Appendix III:	
Digestion, Glucosylated-5-hmC DNA using other GSREs .	12
References .....	13
Contact Information .....	14

## INTRODUCTION

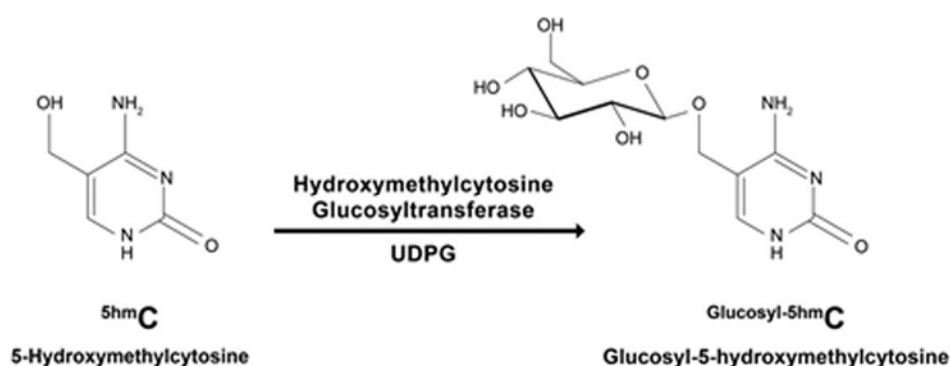
5-Hydroxymethylcytosine (<sup>5</sup>hmC), also known as the “6th base”, is a newly discovered epigenetic modification. Although it was first identified in bacteriophages, its role was thought to be limited to protecting the phage genome from host induced restriction endonucleases<sup>1</sup>. Interestingly, 5-hydroxymethylcytosine has recently been found in embryonic stem cells, the brain, as well as numerous other organs<sup>2,3</sup>. Even though its presence has been confirmed, its biological role remains elusive. It has been proposed that 5-hydroxymethylcytosine may play a role in DNA demethylation (5-methylcytosine) or be involved in another layer of gene expression regulation<sup>3</sup>.

To date, studies of 5-hydroxymethylcytosine has been limited to global quantification or characterization studies. The reason being is that the “standard” for 5-methylcytosine detection, bisulfite DNA analysis, is unable to effectively distinguish 5-methylcytosine from 5-hydroxymethylcytosine.

The **BioArray™ 5-hmC Methylation Kit** allows for sequence specific detection of 5-hydroxymethylcytosine within DNA using a simple and efficient reaction setup.

## ASSAY PRINCIPLE

The highly specific and robust **5-hmC Glucosyltransferase** enzyme, 5-hydroxymethylcytosine in DNA is specifically tagged with a glucose moiety yielding a modified base, glucosyl-5-hydroxymethylcytosine (Figure 1).

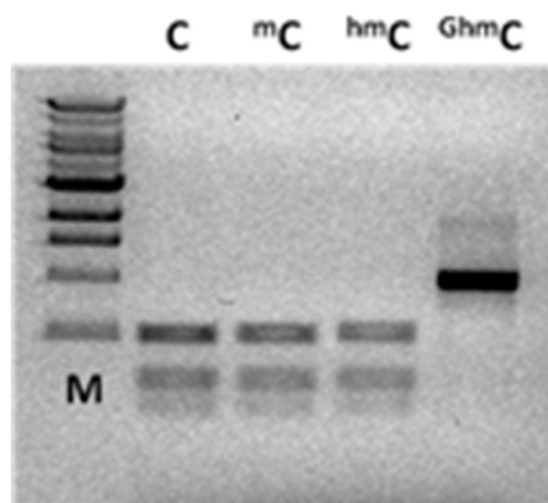


**Figure 1:** 5-hmC Glucosyltransferase transfers a glucose moiety from uridine diphosphoglucose (UDPG) onto preexisting 5-hydroxymethylcytosines within DNA.

After glucosylation of 5-hydroxymethylcytosine, digestion of DNA with “glucosyl-5-hydroxymethylcytosine sensitive” restriction endonucleases (GSREs) allow differentiation of 5-methylcytosine from 5-hydroxymethylcytosine according to the context of a GSRE’s recognition sequence (see **Table 1**). GSREs can efficiently digest DNA when cytosine, 5-methylcytosine, or 5-hydroxymethylcytosine is within their recognition sequence. However, if 5-hydroxymethylcytosine is glucosylated (i.e., glucosyl-5-hydroxymethylcytosine), GSREs can no longer digest the DNA (**Figure 2**). Exploitation of this sensitivity to glucosyl-5-hydroxymethylcytosine facilitates the effective detection 5-hydroxymethylcytosine using a number of downstream applications (e.g., qPCR, NextGen Sequencing, Southern blotting, array, etc.). qPCR is recommended to assess locus-specific context of 5-hydroxymethylcytosine within the DNA Treatment of DNA containing <sup>5hm</sup>C with 5-hmC Glucosyltransferase specifically adds a glucose moiety yielding glucosyl-5-hydroxymethylcytosine. Subsequent digestion with GSREs (see Table 1 below) will specifically cleave DNA having cytosine, 5-methylcytosine, or 5-hydroxymethylcytosine in the recognition sequence. Glucosyl-5-hydroxymethylcytosine in the DNA will remain uncleaved.

GSRE	Recognition Sequence
<b>MspI</b> (included w/ this kit)	C <u><i>CGG</i></u>
<b>GlaI</b>	GCGC
<b>Csp6I</b>	GTAC
<b>HaeIII</b>	GG <u><i>CC</i></u>
<b>Taq<sup>q</sup>I</b>	TCGA
<b>MboI</b>	GATC
<b>McrBC</b>	R <sup>m</sup> C(N <sub>40-3000</sub> )R <sup>m</sup> C

**Table 1: <sup>Glucosyl-5hm</sup>C Sensitive Restriction Endonucleases (GSREs).** Detection of 5-hmC with MspI and HaeIII is only at the indicated “C” position (underlined and italicized). Taq<sup>q</sup>I and MboI display incomplete sensitivity to <sup>Glucosyl-5hm</sup>C. Titration of the enzyme amount and incubation time will be required to yield optimal results.



**Figure 2: Glucosyltransferase (5-hmC) inhibits cleavage of 5hmC DNA by the GSRE Csp6I.** DNA with all cytosines as: unmodified (C), 5-methylcytosine (<sup>m</sup>C), 5-hydroxymethylcytosine (<sup>hm</sup>C) or glucosyl-5-hydroxymethylcytosine (<sup>Ghm</sup>C) was digested with 4 units of Csp6I for 2 hours at 37°C and resolved in a 0.8% w/v agarose/TAE/EtBr gel. “M” is 1kb DNA ladder.

After processing DNA with the BioArray™5-hmC Methylation Kit subsequent analysis of <sup>5hm</sup>C sites can be achieved with qPCR, ultra-deep sequencing, Southern blotting, or microarray.

Additional information on the assay principle and applications can be found in the Appendix I Section of the manual.

## MATERIALS SUPPLIED

BioArray™ 5-hmC Detection Kit	Supply	Storage Temp.
5-hmC Glucosyltransferase (5-hmC)*	100 µL (2 units/µl)	-20 °C
10X 5-hmC Reaction Buffer	1 mL	-20 °C
10X UDPG (Uridine Diphosphoglucose)	600 µL (1mM)	-20 °C
MspI Restriction Enzyme*	150 µL (10 units/µl)	-20 °C
5-hmC Control DNA <sup>(1)</sup>	300 ng	-20 °C
qPCR Primers 1 <sup>(2)</sup>	1 Unit (20 µM )	-20 °C
qPCR Primers 2 <sup>(2)</sup>	1 Unit (20 µM)	-20 °C
5-DNA Clean up Concentrator Kit **	50 preps	RT
<ul style="list-style-type: none"> <li>• DNA Binding</li> <li>• DNA Wash Buffer</li> <li>• DNA Elution Buffer</li> <li>• Spin Column</li> <li>• Collection Tube</li> </ul>		

\*AVOID MULTIPLE FREE THAWING AND LONG TERM STORAGE -80°C.

\*\*ADDITIONAL COMPONENTS PROVIDED

- (1) **5-hmC Control DNA** – A 90bp DNA standard used for qPCR. [10ng/µl] in storage buffer (10mM Tris-HCl; 0.1mM EDTA; pH 8.0) supplied in 30µl volume. See **Appendix** for more details.
- (2) **qPCR Primers 1 and 2** – Control primers specific to 5-hmC Control DNA (above) validated for qPCR. See **Appendix** for detailed information.

## STORAGE

Store reagents suggested conditions. \*RT-Room Temperature

## STABILITY

12 months upon receipt

## PRECAUTIONS

1. Wear appropriate personal protective apparel. Avoid contact with clothes and exposed skin. In case of accidental skin exposure, flush with water immediately.
2. Since some reagents are irritants, wear protective gloves and eye protection. Should be used by trained professionals
3. Interpretation of the results is the sole responsibility of the user.



Reagents require separate storage conditions.

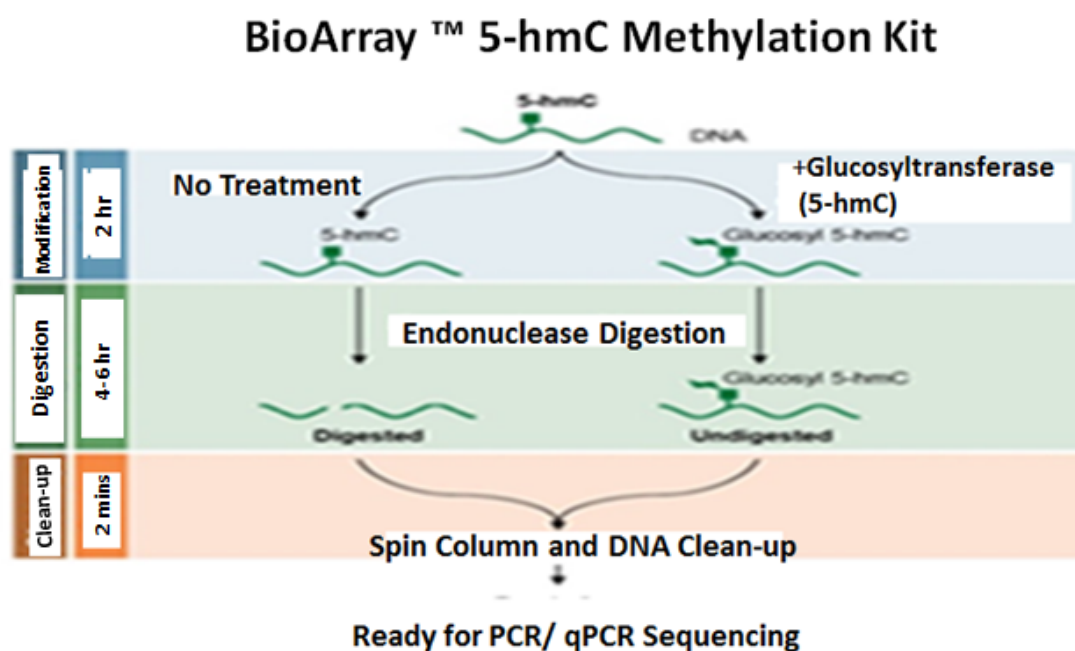
## REAGENT PREPARATION

**NOTE:** Reagents stored at freezing temperatures should be removed prior to experiment and thawed. Reagents should be briefly centrifuged to ensure contents are fully collected.

**5-hmC Control DNA** – A 90bp DNA standard used for qPCR. [10ng/μl] in storage buffer (10mM Tris-HCl; 0.1mM EDTA; Ph 8.0) supplied in 30μl volume. See **Appendix** for more details.

**qPCR Primers 1 and 2** – Control primers specific to **5-hmC Control DNA** (above) validated for qPCR. See **Appendix** for detailed information.

## METHODS AND PROCEDURES



**Figure 3: Overall work flow of the BioArray™ 5-hmC Methylation Kit** (<sup>5hmC</sup> detection in DNA). DNA samples are first modified via glucosylation of <sup>5hmC</sup> within DNA and then digested with MspI. A low elution spin column is featured for clean-up of ultra-pure DNA which is then used in qPCR for locus-specific analysis of <sup>5hmC</sup>.



## PROTOCOL

**Note:** the procedure is also compatible with other GSREs (please see Table 1 and **Appendix**), and other downstream molecular analysis procedures including NextGen sequencing, Southern blotting, array, etc.

### A. Glucosylation Reaction

Divide equal amounts of each DNA to be tested into each of the two separate reaction setups (1 & 2) below. Reaction mixtures containing 5-hmC Control DNA should be set up in parallel (separately) to test DNAs.

#### 1. + Glucosylation Reaction Setup

Test DNA [100-500 ng] or 5-hmC Control DNA [20-50 ng]	x $\mu$ l
10X 5-hmC GT Reaction Buffer	5 $\mu$ l
10X UDPG [1mM]	5 $\mu$ l
5-hmC Glucosyltransferase (2 units/ $\mu$ l)	2 $\mu$ l
ddH <sub>2</sub> O	x $\mu$ l
<b>Total</b>	<b>50 <math>\mu</math>l</b>

**Note:** To ensure the glucosylation reaction is complete, it is recommended that: **a.** A higher enzyme to DNA ratio is used (e.g., use 4 units of 5-hmC Glucosyltransferase:1 $\mu$ g of DNA). **b.** Extend incubation times to  $\geq$ 2 hours at 37°C.

#### 2. Glucosylation Reaction Setup

Test DNA [100-500 ng] or 5-hmC Control DNA [20-50 ng]	x $\mu$ l
10X 5-hmC Reaction Buffer	5 $\mu$ l
10X UDPG [1mM]	5 $\mu$ l
5-hmC GT Enzyme (2 units/ $\mu$ l)	--
ddH <sub>2</sub> O	x $\mu$ l
<b>Total</b>	<b>50 <math>\mu</math>l</b>

Incubate reaction mixtures with DNA at 37°C for  $\geq$  2 hours.

## B. Digestion w/ MspI

**Note:** Alternative GSREs (e.g., Csp6I, HaeIII) can be used in lieu of or together with MspI. Glal is a methylation dependent GSRE. Test DNA *must* be fully methylated with CpG and GpC DNA methyltransferase *in vitro* prior to cleavage with Glal. Glal is compatible with the *single tube reaction format*, but cannot be used in conjunction with other GSREs.

1. Following the incubation from Section **A 1 or 2**, add 30 units of MspI enzyme directly to each (+Glucosylation and -Glucosylation) reaction mixture.
2. Incubate at 37°C for ≥ 4 hours.

**Note:** Extended digestion times and/or additional MspI may be necessary for complete digestion of DNA which is important for subsequent analysis.

## C. Spin Column Clean-up

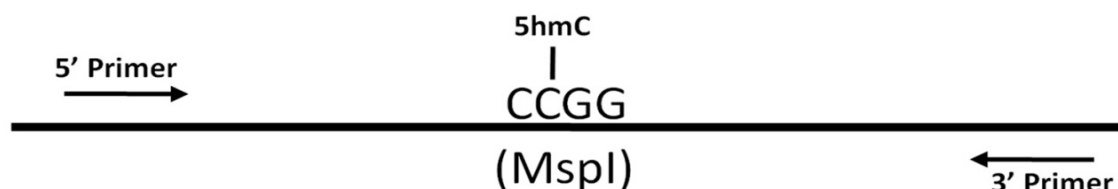
1. Review the components of the **5- DNA Clean up Concentrator Kit** that is included within this kit.
2. Add a 5:1 ratio of **DNA Binding Buffer** to the reaction mixtures (e.g., 250 µl **DNA Binding Buffer** to each 50 µl reaction mixture from Step B)
3. Transfer mixture to the provided **Spin Column** that is placed in the Collection Tube. Centrifuge for 30 seconds (between 10,000-16,000 xg). Discard the flow-through.
4. Add 200µL DNA Wash Buffer to the column. Centrifuge for 30 seconds (between 10,000-16,000 xg). Repeat the wash step.
5. Add ≥6µL DNA Elution Buffer directly to the column matrix and incubate at room temperature for one minute. Transfer the column to a new 1.5mL microcentrifuge tube and centrifuge (between 10,000-16,000 xg) for 30 seconds to elute the DNA.

**Note:** Elution from the column is dependent on pH and temperature. DNA elution Buffer consists of 10mM Tris-HCL pH 8.5, 0.1 mM EDTA. If water is used, make sure that the pH is >6.0. Wait one minute to improve the yield of larger DNA.

6. Following elution of the DNA, qPCR is recommended for locus-specific 5-hmC analysis (see **Step D** below).

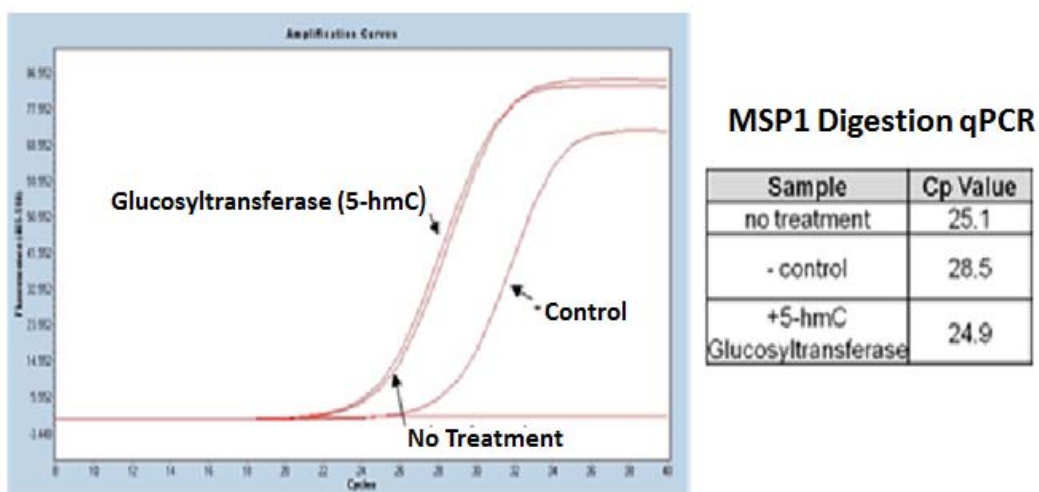
## D. Locus-Specific 5-hmC Identification via qPCR

**Note:** For effective analysis of 5-hmC by qPCR, primers must be designed to flank the GSRE recognition site of interest as shown in **Figure 4**.



**Figure 4:** Design of primers for the area of interest to complete qPCR.

**Note 2:** Each locus requires a proper control for accurate detection of 5-hmC by qPCR as shown in **Figure 5**. This can be achieved by adding test DNA to the qPCR that has not been processed according to **steps A-C** (above).



**Figure 5:** Detection of 5-hmC by qPCR. The “no treatment” DNA will establish the level (i.e., Cp value) that is representative of complete hydroxymethylation at the interrogation site. Conversely, the “- control” (unglucosylated sample) will indicate a level of no hydroxymethylation at the site. The “+Glucosyltransferase (5-hmC)” i.e. glucosylated sample, Cp value should be somewhere between the two, i.e., “no treatment” and “- control”. This will depend on the extent of hydroxymethylation at the interrogation site. *GSRE digested unglucosylated DNA (i.e., “- control”) will always yield a signal albeit at a higher Cp value than the other samples.*

## Appendix I

### 5-hmC Control DNA

The **5-hmC Control DNA** is provided for gauging glucosylation and GSRE reaction efficiencies via qPCR. Incorporated into the **5-hmC Control DNA** are 3 GSRE sites (MspI, HaeIII and Csp6I) which can be used for digestion. The **5-hmC Control DNA** is a 90 bp, double stranded DNA that is partially hemi-5-hydroxymethylated. All cytosine residues in the sequence below that are highlighted in grey (i.e., top strand only) are hydroxymethylated.

#### Sequence of 5-hmC Control DNA:

```

1 ..... 90
caaggatcgctcgcggtctttaGGCCGGtaactgtctgcagctctgagG
TACgcatggattgtaggcgccgacctataccttgtctgcct
gttcttagcgagcgccgagaatCCGGCCattgacagacgtcgagactcC
ATGcgtacctaacatccgcgggcgatgatggaacagacgga
  
```

**Note:** All cytosines in the highlighted sequence are 5-hydroxymethylcytosine. MspI (CCGG), HaeIII (GGCC) and Csp6I (GTAC) sites are capitalized and underlined. MspI and HaeIII sites overlap.

#### Sequence of qPCR Primers:

<b>qPCR Primer 1</b>	5'- caaggatcgctcgcggtcttta -3'
<b>qPCR Primer 2</b>	5'- aggcagacaaggatatagggcg -3'

## Appendix II

### Glucosylation of 5-hmC in DNA

Use the protocol below to completely glucosylate 5-hmC in DNA (without the GSRE digestion and qPCR analysis steps given in the standard protocol of the manual)

**Note:** This can be used for the quantification of global <sup>5hm</sup>C via isotope labeling using Uridine Diphosphate Glucose [Glucose-<sup>14</sup>C(U)].

### Glucosylation Reaction Setup

DNA [100-500 ng]	10 µl
10X 5-hmC GT Reaction Buffer	5 µl
10X UDPG [1mM]	5 µl
5-hmC Glucosyltransferase Enzyme (2 units/µl)	2 µl
ddH <sub>2</sub> O	28 µl
<b>Total</b>	<b>50 µl</b>

5-hmC in DNA should be fully glucosylated after incubating the reaction at 37°C for ≥ 2 hours. DNA can be recovered using the clean-up procedure in the protocol (**Step C in Methods and Procedures**)

## Appendix III

### Digestion of Glucosylated-5-hmC DNA using other GSREs

For the GSREs MspI, Glal, Csp6I, and HaeIII, glucosylation followed by digestion can be performed (i.e. single tube reaction format). However, TaqI, MboI, and McrBC are incompatible with the *single tube reaction format* and the protocol given below should be used with these enzymes instead.

#### Step 1 – Glucosylation Reaction

Divide equal amounts of each DNA to be tested into each of the two separate reaction setups (**A & B**) below. Reaction mixtures containing **5-hmC Control DNA** should be set up in parallel (separately) to test DNAs.

##### **A. + Glucosylation Reaction Setup**

Test DNA [100-500 ng] or 5-hmC Control DNA [20-50 ng]	x $\mu$ l
10X 5-hmC Reaction Buffer	5 $\mu$ l
10X UDPG [1mM]	5 $\mu$ l
5-hmC Glucosyltransferase (5-hmC ) (2 units/ $\mu$ l)	2 $\mu$ l
ddH <sub>2</sub> O	x $\mu$ l
<b>Total</b>	<b>50 <math>\mu</math>l</b>

##### **B. Glucosylation Reaction Setup**

Test DNA [100-500 ng] or 5-hmC Control DNA [20-50 ng]	x $\mu$ l
10X 5-hmC Reaction Buffer	5 $\mu$ l
10X UDPG [1mM]	5 $\mu$ l
5-hmC Glucosyltransferase (5-hmC ) (2 units/ $\mu$ l)	--
ddH <sub>2</sub> O	x $\mu$ l
<b>Total</b>	<b>50 <math>\mu</math>l</b>

Incubate reaction mixtures with DNA at 37°C for  $\geq$  2 hours.

## Step 2 – Spin Column Clean-up

- a. Add a 5:1 ratio of **DNA Binding Buffer** to the reaction mixtures (e.g., 250µl **DNA Binding Buffer** to each 50µl reaction mixture).
- b. Transfer mixture to the provided **Spin Column** that is placed in the Collection Tube. Centrifuge for 30 seconds (between 10,000-16,000 xg). Discard the flow-through.
- c. Add 200µL DNA Wash Buffer to the column. Centrifuge for 30 seconds (between 10,000-16,000 xg). Repeat the wash step.
- d. Add ≥6µL DNA Elution Buffer directly to the column matrix and incubate at room temperature for one minute. Transfer the column to a new 1.5mL microcentrifuge tube and centrifuge (between 10,000-16,000 xg) for 30 seconds to elute the DNA.

## Step 3 – GSRE Digestion

Eluted DNA from the previous step should be digested with GSRE following supplier's recommended conditions.

**Note:** *Over digestion of DNA is recommended. Extended digestion times and/or additional units of GSRE may be necessary for complete digestion of DNA which is important for subsequent analysis.*

## Step 4 – Spin Column Clean-up

Repeat Step 2 from this Appendix Section. Following elution of the DNA, qPCR is recommended for locus-specific 5-hmC analysis.

## REFERENCES

1. Wyatt GR, Cohen SS (1952) . A new pyrimidine base from bacteriophage nucleic acids. Nature 170 (4338) 1072-3, 1952.
2. Riaucionis S., Heintz N., (2009). The nuclear DNA base 5-hydroxymethylcytosine is present in Purkinje neurons and the brain. Science 324 (5929): 929-30.
3. Tahiliani M., et al. (2009). "Conversion of 5- methylcytosine to 5- hydroxymethylcytosine in mammalian DNA by MLL partner TET1". Science 324 (5929): 930–35
4. Huang Y., et al. (2010). "The behaviour of 5-hydroxymethylcytosine in bisulfite sequencing". PLoS One. 2010 Jan 26;5(1)
5. Szwagierczak A., et al. (2010), "Sensitive enzymatic quantification of 5- hydroxymethylcytosine in genomic DNA" Nucleic Acids Res.



# Product Manual

## **GLOBAL HEADQUARTERS**

Enzo Life Sciences Inc.  
10 Executive Boulevard  
Farmingdale, NY 11735  
Toll-Free: 1.800.942.0430  
Phone: 631.694.7070  
Fax: 631.694.7501  
info-usa@enzolifesciences.com

## **EUROPE/ASIA**

Enzo Life Sciences (ELS) AG  
Industriestrasse 17  
CH-4415 Lausen  
Switzerland  
Phone: +41/0 61 926 89 89  
Fax: +41/0 61 926 89 79  
info-ch@enzolifesciences.com

For local distributors and detailed product information visit us online:  
[www.enzolifesciences.com](http://www.enzolifesciences.com)