

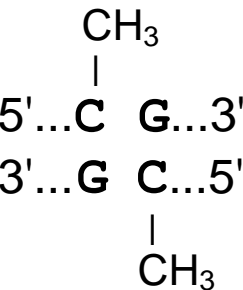


PRODUCT INFORMATION

CpG Methyltransferase (M.SssI)

Pub. No. MAN0011994
Rev. Date 15.July.2016 (Rev. B.00)

#_ _
Lot _ _ Expiry Date _ _



Components	#EM0821
M.SssI	25 µL
10X M.SssI Buffer	1 mL
50X (5 mM) SAM	0.1 mL

Store at -20 °C

www.thermofisher.com

For Research Use Only. Not for use in diagnostic procedures.

Description

CpG Methyltransferase (M.SssI) is one of the basic tools used in epigenetic studies. The enzyme methylates the C⁵ position on the base moiety of all cytosine nucleotides contained in unmethylated or hemimethylated double stranded DNA in a 5'-CpG-3' context. Thermo Scientific M.SssI is specifically formulated for fast reaction times without compromising the reaction efficiency. The enzyme completes modification of all CpGs in 15 min at 37 °C. In addition, the enzyme has been specifically validated for use with genomic DNA – the primary substrate in epigenetic studies. M.SssI is supplied with an optimized 10X reaction buffer and 50X S-adenosylmethionine (SAM) as a cofactor.

Storage Buffer

CpG Methyltransferase (M.SssI) is supplied in: 10 mM potassium phosphate (pH 7.0 at 25 °C), 400 mM KCl, 1 mM DTT, 1 mM EDTA, 0.2 mg/mL BSA and 50% (v/v) glycerol.

Features

- **High efficiency** - complete *in vitro* methylation of all CpG sequences in non-methylated and hemi-methylated DNA.
- **Fast** - reaction is completed in 15 min at 37 °C.
- **Stable** - at -20 °C for 2 years.

Applications

- Epigenetics studies.
- *In vitro* methylation of DNA for methylation analysis.
- Inhibition of endonucleases with overlapping CpG sequence recognition.
- [³H]-labeling of DNA.

Enzyme formulation

One μ L of M.SssI protects 1 μ g of genomic DNA from digestion with HpaII in 15 min at 37 °C in 50 μ L of recommended reaction buffer.

CERTIFICATE OF ANALYSIS

Prolonged incubation

No detectable change in DNA fragment band pattern observed following agarose gel electrophoresis of SmaI linearized pUC19 DNA incubated for 16 hours with 1 μ L of M.SssI.

Labeled Oligonucleotide (LO) Assay

No detectable degradation after incubation of single-stranded or double-stranded radiolabeled oligonucleotides with M.SssI.

Quality authorized by:

 Jurgita Zilinskiene

Protocol

- Assemble the following reaction at room temperature:

Nuclease-free water	to 20 μ L
10X M.SssI Buffer	2 μ L
50X SAM	0.4 μ L
DNA	up to 1 μ g
M.SssI	1 μ L
Total volume	20 μ L

- Mix gently and spin down for a few seconds.
- Incubate at 37 °C for 15 min.
- Stop the reaction by heating at 65 °C for 20 min.

DNA can be purified by phenol extraction followed by ethanol precipitation or by using a spin column kit.

LIMITED USE LABEL LICENSE: Internal Research and Development Use Only.

The purchase of this product conveys to the buyer the limited, non-exclusive, non-transferable right (without the right to resell, repackage, or further sublicense) to use this product for internal research and development purposes. No other license is granted to the buyer whether expressly, by implication, by estoppel or otherwise. In particular, the purchase of the product does not include or carry any right or license to use, develop, or otherwise exploit this product commercially and no rights are conveyed to the buyer to use the product or components of the product for purposes including but not limited to provision of services to a third party, generation of commercial databases or clinical diagnostics. This product is sold pursuant to authorization from Thermo Fisher Scientific and Thermo Fisher Scientific reserves all other rights. For information on purchasing a license for uses other than internal research and development purposes, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies Inc., 5791 Van Allen Way, Carlsbad, California 92008.

PRODUCT USE LIMITATION

This product is developed, designed and sold exclusively *for research purposes and in vitro use only*. The product was not tested for use in diagnostics or for drug development, nor is it suitable for administration to humans or animals.

Please refer to www.thermofisher.com for Material Safety Data Sheet of the product.

© 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.