



Innovating Epigenetics Solutions

# Premium WGBS Indexes

## For Whole Genome Bisulfite Sequencing

**Cat. No.** C05010032 (48 rxns, 6 indexes)  
C05010033 (96 rxns, 12 indexes)  
C05010034 (192 rxns, 24 indexes)



Please read this manual carefully  
before starting your experiment

# Contents

---

Introduction	4
Kit materials	5
Remarks before starting	5
Oligonucleotide sequences	6

# Introduction

---

The **Premium Whole Genome Bisulfite Sequencing (WGBS) Indexes** have been designed to prepare multiplexed single and paired-end bisulfite converted DNA libraries for sequencing using Illumina® platforms. The index and flow cell binding sequence are designed within the Premium WGBS adapters and are attached onto the sample insert during adapter ligation. Pooling with Premium WGBS Indexes allows for multiplexing several samples in a single flow cell.

The Premium WGBS Indexes are compatible with Diagenode's Premium WGBS kit (Cat. No. C02030034) and suitable for performing Whole Genome Bisulfite Sequencing or ChIP-Bisulfite-Sequencing.

# Kit materials

---

The Premium WGBS Indexes are available in three different formats containing 6, 12 or 24 Methylated Barcoded Adapters with enough material for 8 reactions each. The shelf life of each reagent is 12 months when stored at -20°C.

Caution: Store the components at the indicated temperature upon receipt.

Table 1a. Components supplied with the Premium WGBS

Description	Quantity	Storage
Premium WGBS Indexes (25µM): 6, 12 or 24 tubes (purple)	1.5 mL	-20°C
Premium WGBS Primer Mix (green)	56 µL	-20°C

Caution: The Premium WGBS Barcoded Adapters are supplied in duplex form. Do not heat the adapters above room temperature.

## Remarks before starting

---

- It is highly recommended to use the supplied Premium WGBS Primer Mix during PCR amplification. Inadvertent use of an incorrect primer sequence can potentially result in elimination of the index.
- Do not heat the Premium WGBS Indexes above room temperature.

# Oligonucleotide sequences

PCR Primer 1	5'AATGATACGGCGACCACCGAGATCTACAC
PCR Primer 2	5'CAAGCAGAAGACGGCATAACGAGAT
Bisulfite-seq Adapters	5'AATGATACGGCGACCACCGAGATCTACACTCTTTCCCTACACGACGCTCTTCCGATCT 5'GATCGGAAGAGCACACGTCTGAACTCCAGTCACXXXXX <sup>1</sup> ATCTCGTATGCCGTCTTCTGCTTG

<sup>1</sup>XXXXXX denotes the index region of the adapter. The index sequences contained in each adapter are listed below.

Premium WGBS Indexes - 6	
Premium WGBS Index 24	GGTAGC
Premium WGBS Index 26	ATGAGC
Premium WGBS Index 31	CACGAT
Premium WGBS Index 38	CTAGCT
Premium WGBS Index 41	GCGCTA
Premium WGBS Index 44	TATAAT

Premium WGBS Indexes - 12	
Premium WGBS Index 24	GGTAGC
Premium WGBS Index 26	ATGAGC
Premium WGBS Index 28	CAAAAG
Premium WGBS Index 29	CAACTA
Premium WGBS Index 30	CACCGG
Premium WGBS Index 31	CACGAT
Premium WGBS Index 32	CACTCA
Premium WGBS Index 33	CAGGCG
Premium WGBS Index 34	CATGGC
Premium WGBS Index 38	CTAGCT
Premium WGBS Index 41	GCGCTA
Premium WGBS Index 44	TATAAT

Premium WGBS Indexes - 24	
Premium WGBS Index 17	GTAGAG
Premium WGBS Index 24	GGTAGC
Premium WGBS Index 26	ATGAGC
Premium WGBS Index 28	CAAAAG
Premium WGBS Index 29	CAACTA
Premium WGBS Index 30	CACCGG
Premium WGBS Index 31	CACGAT
Premium WGBS Index 32	CACTCA
Premium WGBS Index 33	CAGGCG
Premium WGBS Index 34	CATGGC
Premium WGBS Index 35	CATTTT
Premium WGBS Index 36	CCAACA
Premium WGBS Index 37	CGGAAT
Premium WGBS Index 38	CTAGCT
Premium WGBS Index 39	CTATAC
Premium WGBS Index 40	CTCAGA
Premium WGBS Index 41	GCGCTA
Premium WGBS Index 42	TAATCG
Premium WGBS Index 43	TACAGC
Premium WGBS Index 44	TATAAT
Premium WGBS Index 45	TCATTC
Premium WGBS Index 46	TCCCGA
Premium WGBS Index 47	TCGAAG
Premium WGBS Index 48	TCGGCA

Caution: For low level multiplexing use the following barcodes:

For 2 barcodes: 24 and 31 or 38 and 41

For 3 barcodes: 24, 31 and any other barcode or 38, 41 and any other barcode

## **FOR RESEARCH USE ONLY.**

### **Not intended for any animal or human therapeutic or diagnostic use.**

© 2017 Diagenode SA. All rights reserved. No part of this publication may be reproduced, transmitted, transcribed, stored in retrieval systems, or translated into any language or computer language, in any form or by any means: electronic, mechanical, magnetic, optical, chemical, manual, or otherwise, without prior written permission from Diagenode SA (hereinafter, "Diagenode"). The information in this guide is subject to change without notice. Diagenode and/or its affiliates reserve the right to change products and services at any time to incorporate the latest technological developments. Although this guide has been prepared with every precaution to ensure accuracy, Diagenode and/or its affiliates assume no liability for any errors or omissions, nor for any damages resulting from the application or use of this information. Diagenode welcomes customer input on corrections and suggestions for improvement.

### **NOTICE TO PURCHASER LIMITED LICENSE**

The information provided herein is owned by Diagenode and/or its affiliates. Subject to the terms and conditions that govern your use of such products and information, Diagenode and/or its affiliates grant you a nonexclusive, nontransferable, non-sublicensable license to use such products and information only in accordance with the manuals and written instructions provided by Diagenode and/or its affiliates. You understand and agree that except as expressly set forth in the terms and conditions governing your use of such products, that no right or license to any patent or other intellectual property owned or licensable by Diagenode and/or its affiliates is conveyed or implied by providing these products. In particular, no right or license is conveyed or implied to use these products in combination with any product not provided or licensed to you by Diagenode and/or its affiliates for such use. Limited Use Label License: Research Use Only The purchase of this product conveys to the purchaser the limited, non-transferable right to use the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial applications of any kind, including, without limitation, quality control and commercial services such as reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact [info@diagenode.com](mailto:info@diagenode.com).

### **TRADEMARKS**

The trademarks mentioned herein are the property of Diagenode or their respective owners. Bioanalyzer is a trademark of Agilent Technologies, Inc. Agencourt and AMPure® are registered trademarks of Beckman Coulter, Inc. Illumina® is a registered trademark of Illumina® Inc; Qubit is a registered trademark of Life Technologies Corporation.

[www.diagenode.com](http://www.diagenode.com)