

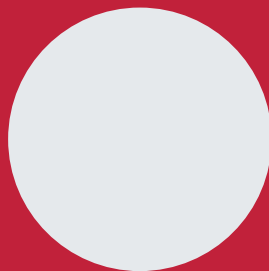
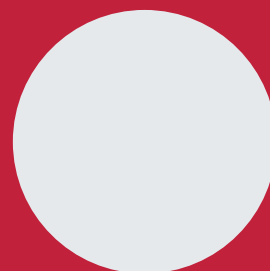


Innovating Epigenetic Solutions

Premium WGBS Indexes

For Whole Genome Bisulfite Sequencing

Cat. No. **C05010032** (48 rxns, 6 indices)
C05010032 (96 rxns, 12 indices)
C05010032 (192 rxns, 24 indices)



Content

Introduction	4
Kit materials	5
Remarks before starting	6
Oligonucleotide sequences	11

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 Barcodes kit

Description	Quantity	Storage
WGBS Adapters (25µM): 6, 12 or 24 tubes (purple)	20 µl	-20°C
WGBS Primer Mix (green)t	96 µg	-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 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 WGBS Indexes above room temperature.

Oligonucleotide sequences

PCR Primer 1	5'AATGATACGGCGACCACCGAGATCTACAC
PCR Primer 2	5'CAAGCAGAAGACGGCATACGAGAT
Bisulfite-seq Adapters	5'AATGATACGGCGACCACCGAGATCTACACTCTTTCCCTACACGACGCTCTTCCGATCT 5'GATCGGAAGAGCACACGTCTGAACTCCAGTCACXXXXX1ATCTCGTATGCCGTCTTCTGCTTG

1 XXXXX denotes the index region of the adapter. The index sequences contained in each adapter are listed below.

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

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

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

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.**

© 2016 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.

TRADEMARKS

The trademarks mentioned herein are the property of Diagenode or their respective owners. Illumina is a registered trademark of Illumina Inc.



Innovating Epigenetic Solutions

**DIAGENODE
HEADQUARTERS**

**DIAGENODE S.A.
BELGIUM | EUROPE**

LIEGE SCIENCE PARK
Rue Bois Saint-Jean, 3
4102 Seraing - Belgium
Tel: +32 4 364 20 50
Fax: +32 4 364 20 51
orders@diagenode.com
info@diagenode.com

**DIAGENODE INC.
USA | NORTH AMERICA**

400 Morris Avenue, Suite #101
Denville, NJ 07834
Tel: +1 862 209-4680
Fax: +1 862 209-4681
orders.na@diagenode.com
info.na@diagenode.com

**FOR A COMPLETE LISTING OF
DIAGENODE'S INTERNATIONAL
DISTRIBUTORS VISIT:**

[http://www.diagenode.com/company/
distributors.php](http://www.diagenode.com/company/distributors.php)
For rest of the world, please contact
Diagenode sa.
www.diagenode.com