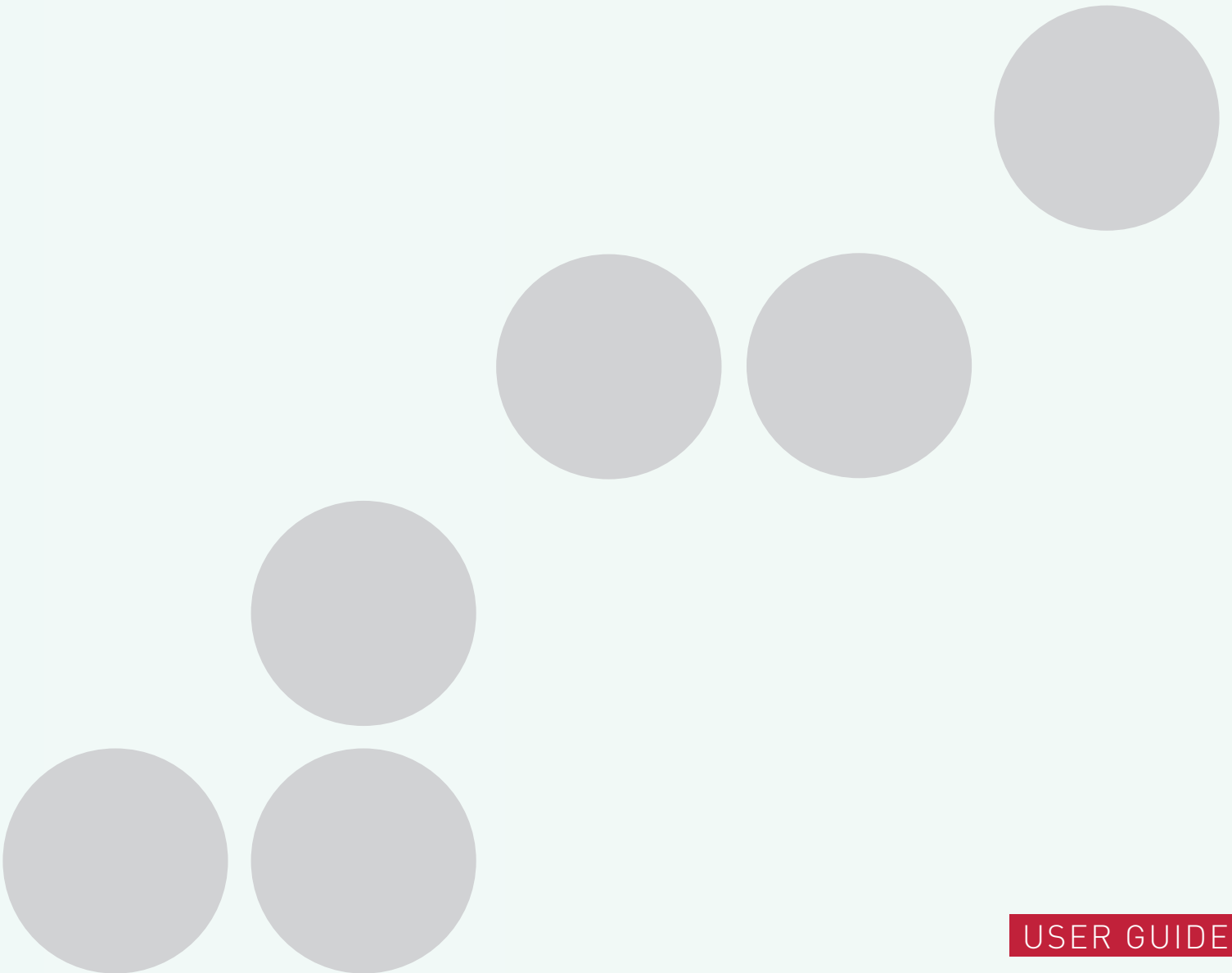




Innovating Epigenetics Solutions

24 SI for ChIPmentation

Cat. No. C01011031 (24 rxns)





Please read this manual carefully
before starting your experiment

Contents

Introduction	4
Indexes	5
Multiplexing and index pooling	6
Related products	6

Introduction

The **24 SI for ChIPmentation** includes 24 single primer indexes allowing the multiplexing of up to 24 samples for sequencing on Illumina platforms. This set of indexes is designed to be used with TAG Kit for ChIPmentation (Diagenode, Cat. No. C01011030), but it is also compatible with other tagmentation-based library preparation protocols, such as ATAC-Seq or CUT&Tag technologies. For ChIPmentation protocol please follow the manual of the TAG Kit for ChIPmentation.

Components supplied with the 24 SI for ChIPmentation

NOTE: Upon receipt, store the components at the indicated temperatures.

Table 1. Components stored at -20°C

Component	Cap color	Volume	Storage
Universal primer	transparent	44 µl	-20°C
Primer Index 1	transparent	12 µl	-20°C
Primer Index 2	none	4,5 µl	-20°C
Primer Index 3	none	4,5 µl	-20°C
Primer Index 4	none	4,5 µl	-20°C
Primer Index 5	none	4,5 µl	-20°C
Primer Index 6	none	4,5 µl	-20°C
Primer Index 7	none	4,5 µl	-20°C
Primer Index 8	none	4,5 µl	-20°C
Primer Index 9	none	4,5 µl	-20°C
Primer Index 10	none	4,5 µl	-20°C
Primer Index 11	none	4,5 µl	-20°C
Primer Index 12	none	4,5 µl	-20°C
Primer Index 13	none	4,5 µl	-20°C
Primer Index 14	none	4,5 µl	-20°C
Primer Index 15	none	4,5 µl	-20°C
Primer Index 16	none	4,5 µl	-20°C
Primer Index 17	none	4,5 µl	-20°C
Primer Index 18	none	4,5 µl	-20°C
Primer Index 19	none	4,5 µl	-20°C
Primer Index 20	none	4,5 µl	-20°C
Primer Index 21	none	4,5 µl	-20°C
Primer Index 22	none	4,5 µl	-20°C
Primer Index 23	none	4,5 µl	-20°C
Primer Index 24	none	4,5 µl	-20°C

Indexes

The primer indexes for ChIPmentation contains 8-nucleotide long i7 indexes. Each primer index is provided with enough volume for 3 reactions.

Table 2. List of the 24 primer indexes with their indexes sequences

Reagent	Index Sequence
Primer Index 1	TAAGGCGA
Primer Index 2	CGTACTAG
Primer Index 3	AGGCAGAA
Primer Index 4	TCCTGAGC
Primer Index 5	GGACTCCT
Primer Index 6	TAGGCATG
Primer Index 7	CTCTCTAC
Primer Index 8	CAGAGAGG
Primer Index 9	GCTACGCT
Primer Index 10	CGAGGCTG
Primer Index 11	AAGAGGCA
Primer Index 12	GTAGAGGA
Primer Index 13	GTCGTGAT
Primer Index 14	ACCACTGT
Primer Index 15	TGGATCTG
Primer Index 16	CCGTTTGT
Primer Index 17	TGCTGGGT
Primer Index 18	GAGGGGTT
Primer Index 19	AGGTTGGG
Primer Index 20	GTGTGGTG
Primer Index 21	TGGTTTTC
Primer Index 22	TGGTCACA
Primer Index 23	TTGACCCT
Primer Index 24	CCACTCCT

Multiplexing and index pooling

When multiplexing less than 5 samples together the choice of the index is important to ensure a good quality sequencing. We recommend the following combinations when low multiplexing of libraries is required:

Multiplexing level	Index combination
2 samples	Index 1 - Index 2 Index 2 - Index 4
3 samples	Index 1 - Index 2 - Index 4 Index 3 - Index 5 - Index 6 Index 10 - Index 11 - Index 12 Index 13 - Index 14 - Index 22 2-plex option with any other index
4 samples	Index 8 - Index 9 - Index 11 - Index 12 Index 8 - Index 12 - Index 21 - Index 24 Index 1 - Index 2 - Index 3 - Index 4 Index 4 - Index 7 - Index 19 - Index 23 3-plex option with any other index

For 5 - 11-plex pools use 4-plex options with any other available indexes. If a higher multiplexing degree is required (> 12), any combination is possible regardless of the index chosen.

Related products

Product	Cat. No.
TAG Kit for ChIPmentation	C01011030
iDeal ChIP-seq Kit for Histones	C01010051
iDeal ChIP-seq Kit for Transcription Factors	C01010055
True MicroChIP Kit	C01010130
DiaMag 0.2ml magnetic rack	B04000001
DiaMag 1.5ml magnetic rack	B04000003
DiaMag Rotator	B05000001
Bioruptor Pico®	B01060001
IP-Star® Compact Automated System	B03000002

FOR RESEARCH USE ONLY.

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

© 2019 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. 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