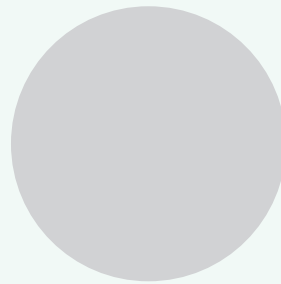
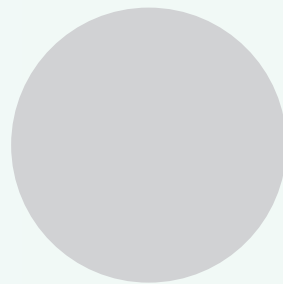
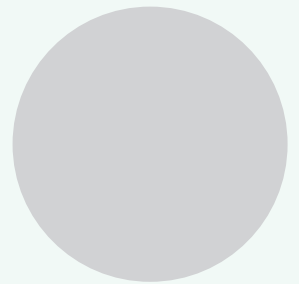




Innovating Epigenetics Solutions

# Primer indexes for tagmented libraries

- Cat. No.** C01011033 – 8 SI for tagmented libraries  
C01011032 – 24 SI for tagmented libraries  
C01011035 – 8 UDI for tagmented libraries  
C01011034 – 24 UDI for tagmented libraries – Set I  
C01011036 – 24 UDI for tagmented libraries – Set II





Please read this manual carefully  
before starting your experiment

# Contents

---

<b>Introduction</b>	<b>5</b>
<b>8 SI for tagmented libraries</b> (Cat. No. C0101133)	<b>6</b>
Multiplexing and index pooling	7
<b>24 SI for tagmented libraries</b> (Cat. No. C01011032)	<b>8</b>
Multiplexing and index pooling	10
<b>8 UDI for tagmented libraries</b> (Cat. No. C01011035)	<b>11</b>
Multiplexing and index pooling	12
<b>24 UDI for tagmented libraries – Set I</b> (Cat. No. C01011034)	<b>13</b>
Multiplexing and index pooling	15
<b>24 UDI for tagmented libraries – Set II</b> (Cat. No. C01011036)	<b>16</b>
<b>Related products</b>	<b>18</b>



# Introduction

---

The primer indexes for tagmented libraries are PCR primers targeting the Nextera sequencing adaptors, previously incorporated in the libraries by tagmentation. The structure of final library is shown in figure 1.

---

## A – Single index libraries:

5'AATGATACGGCGACCACCGAGATCTACACTCGTCGGCAGCGTCAGATGTGTATAAGAGACAG – insert – CTGTCTCTTATACACATCTCCGAGCCCACGAGAC[i7index]ATCTCGTATGCCGTCTTCTGCTTG 3'

## B – Dual index libraries:

5' AATGATACGGCGACCACCGAGATCTACAC[i5 index]TCGTCTCGGCAGCGTCAGATGTGTATAAGAGACAG – insert – CTGTCTCTTATACACATCTCCGAGCCCACGAGAC[i7 index]ATCTCGTATGCCGTCTTCTGCTTG 3'

---

**Figure 1.** Structure of tagmented libraries after final amplification with Diagenode primer indexes. A) Single-index libraries. B) Dual-index libraries.

Each primer index is bearing a unique index (in red in figure 1) in order to identify each library before pooling them for sequencing in the same lane. To give flexibility in the experimental design and answer different needs several products are available:

- **8 SI for tagmented libraries** (Cat. No. C01011033)
- **24 SI for tagmented libraries** (Cat. No. C01011032)
- **8 UDI for tagmented libraries** (Cat. No. C01011035)
- **24 UDI for tagmented libraries – Set I** (Cat. No. C01011034)
- **24 UDI for tagmented libraries – Set II** (Cat. No. C01011036)

They are all compatible with any Nextera-based libraries such as the one generated with ChIPmentation, ATAC-seq or CUT&Tag technologies.

# 8 SI for tagmented libraries

Cat. No. C01011033

The 8 SI for tagmented libraries includes 8 primer pairs for single indexing allowing the multiplexing of up to 8 samples for sequencing on Illumina platforms. This set of indexes is designed to be used with  $\mu$ ChIPmentation Kit for Histones (Cat. No. C01011012) but it is also compatible with other tagmentation-based library preparation protocols, such as ATAC-seq or CUT&Tag technologies.

The primer indexes for tagmented libraries contain 8-nucleotide long i7 indexes.

Table 1. Components supplied with the kit 8 SI for tagmented libraries with their indexes sequences

Component	i7 Index Sequence	Cap color	Volume	Storage
Primer pair SI 1	TAAGGCGA	transparent	24 $\mu$ l	-20°C
Primer pair SI 2	CGTACTAG	none	9 $\mu$ l	-20°C
Primer pair SI 3	AGGCAGAA	none	9 $\mu$ l	-20°C
Primer pair SI 4	TCCTGAGC	none	9 $\mu$ l	-20°C
Primer pair SI 5	GGACTCCT	none	9 $\mu$ l	-20°C
Primer pair SI 6	TAGGCATG	none	9 $\mu$ l	-20°C
Primer pair SI 7	CTCTCTAC	none	9 $\mu$ l	-20°C
Primer pair SI 8	CAGAGAGG	none	9 $\mu$ l	-20°C

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

# 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 SI 1 - Index SI 2 Index SI 2 - Index SI 4
3 samples	Index SI 1 - Index SI 2 - Index SI 4 Index SI 3 - Index SI 5 - Index SI 6 2-plex option with any other index
4 samples	Index SI 1 - Index SI 2 - Index SI 3 - Index SI 4 3-plex option with any other index

For 5 - 8-plex pools use 4-plex options with any other available indexes.

# 24 SI for tagmented libraries

Cat. No. C01011032

The 24 SI for tagmented libraries includes 24 primer pairs for single-indexing allowing the multiplexing of up to 24 samples for sequencing on Illumina platforms. This set of indexes is designed to be used with  $\mu$ ChIPmentation for Histones (Cat. No. C01011011) but it is also compatible with other tagmentation-based library preparation protocols, such as ATAC-Seq or CUT&Tag technologies.

The primer indexes for tagmented libraries contain 8-nucleotide long i7 indexes.

**Table 1.** Components supplied with the 24 SI for tagmented libraries and their indexes sequences

Component	i7 Index Sequence	Cap color	Volume	Storage
Primer pair SI 1	TAAGGCGA	transparent	24 $\mu$ l	-20°C
Primer pair SI 2	CGTACTAG	none	9 $\mu$ l	-20°C
Primer pair SI 3	AGGCAGAA	none	9 $\mu$ l	-20°C
Primer pair SI 4	TCCTGAGC	none	9 $\mu$ l	-20°C
Primer pair SI 5	GGACTCCT	none	9 $\mu$ l	-20°C
Primer pair SI 6	TAGGCATG	none	9 $\mu$ l	-20°C
Primer pair SI 7	CTCTCTAC	none	9 $\mu$ l	-20°C
Primer pair SI 8	CAGAGAGG	none	9 $\mu$ l	-20°C
Primer pair SI 9	GCTACGCT	none	9 $\mu$ l	-20°C
Primer pair SI 10	CGAGGCTG	none	9 $\mu$ l	-20°C
Primer pair SI 11	AAGAGGCA	none	9 $\mu$ l	-20°C
Primer pair SI 12	GTAGAGGA	none	9 $\mu$ l	-20°C
Primer pair SI 13	GTCGTGAT	none	9 $\mu$ l	-20°C
Primer pair SI 14	ACCACTGT	none	9 $\mu$ l	-20°C
Primer pair SI 15	TGGATCTG	none	9 $\mu$ l	-20°C
Primer pair SI 16	CCGTTTGT	none	9 $\mu$ l	-20°C
Primer pair SI 17	TGCTGGGT	none	9 $\mu$ l	-20°C
Primer pair SI 18	GAGGGGTT	none	9 $\mu$ l	-20°C
Primer pair SI 19	AGGTTGGG	none	9 $\mu$ l	-20°C



Component	i7 Index Sequence	Cap color	Volume	Storage
Primer pair SI 20	GTGTGGTG	none	9 $\mu$ l	-20°C
Primer pair SI 21	TGGGTTTC	none	9 $\mu$ l	-20°C
Primer pair SI 22	TGGTCACA	none	9 $\mu$ l	-20°C
Primer pair SI 23	TTGACCCT	none	9 $\mu$ l	-20°C
Primer pair SI 24	CCACTCCT	none	9 $\mu$ l	-20°C

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

# 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 SI 1 – Index SI 2 Index SI 2 - Index SI 4
3 samples	Index SI 1 - Index SI 2 - Index SI 4 Index SI 3 - Index SI 5 - Index SI 6 Index SI 10 - Index SI 11 - Index SI 12 Index SI 13 - Index SI 14 - Index SI 22 2-plex option with any other index
4 samples	Index SI 8 – Index SI 9 – Index SI 11 – Index SI 12 Index SI 8 - Index SI 12 – Index SI 21 – Index SI 24 Index SI 1 – Index SI 2 - Index SI 3 – Index SI 4 Index SI 4 - Index SI 7 - Index SI 19 - Index SI 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.

# 8 UDI for tagmented libraries

Cat. No. C01011035

The 8 UDI (Unique dual indexes) for tagmented libraries provides combinations of barcodes where each barcode is uniquely attributed to one sample. This is a great tool to identify mistakes during index sequencing. Indeed, a phenomenon, known as index hopping, can lead to misattribution of some reads to the wrong sample. This is particularly frequent with the NovaSeq6000, and thus the use of Unique Dual-Indexing (UDI) is therefore highly recommended when using this sequencer.

The 8 UDI for tagmented libraries includes 8 primer pairs for unique dual-indexing allowing the multiplexing of up to 8 samples for sequencing on Illumina platforms. This set of indexes is designed to be used with  $\mu$ ChIPmentation Kit for Histones (Cat. No. C01011012) but it is also compatible with other tagmentation-based library preparation protocols, such as ATAC-seq or CUT&Tag technologies.

The primer indexes for tagmented libraries contain 8-nucleotide long i7 and i5 indexes.

**Table 1.** Components supplied with the kit 8 UDI for tagmented libraries with their indexes sequences

Component	i7 Index Sequence	i5 Index Sequence	Cap color	Volume	Storage
Primer pair UDI 1	AACCGCGG	AGCGCTAG	transparent	12 $\mu$ l	-20°C
Primer pair UDI 2	GGTTATAA	GATATCGA	none	5 $\mu$ l	-20°C
Primer pair UDI 3	CCAAGTCC	CGCAGACG	none	5 $\mu$ l	-20°C
Primer pair UDI 4	TTGGACTT	TATGAGTA	none	5 $\mu$ l	-20°C
Primer pair UDI 5	CAGTGGAT	AGGTGCGT	none	5 $\mu$ l	-20°C
Primer pair UDI 6	TGACAAGC	GAACATAC	none	5 $\mu$ l	-20°C
Primer pair UDI 7	CTAGCTTG	ACATAGCG	none	5 $\mu$ l	-20°C
Primer pair UDI 8	TCGATCCA	GTGCGATA	none	5 $\mu$ l	-20°C

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

# 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	UDI 1 - UDI 2 UDI 1 - UDI 8 UDI 2 - UDI 7 UDI 3 - UDI 4 UDI 5 - UDI 6 UDI 7 - UDI 8
3 samples	UDI 1 - UDI 2 - UDI 7 UDI 1 - UDI 2 - UDI 8 UDI 1 - UDI 7 - UDI 8 UDI 2 - UDI 7 - UDI 8 UDI 5 - UDI 6 - UDI 7 UDI 5 - UDI 6 - UDI 8 UDI 5 - UDI 7 - UDI 8 UDI 6 - UDI 7 - UDI 8
4 samples	3-plex option with any other index

For 5 - 8-plex pools use 4-plex options with any other available indexes.

# 24 UDI for tagmented libraries - Set I

---

Cat. No. C01011034

The 24 UDI (Unique dual indexes) for tagmented libraries – Set I provides combinations of barcodes where each barcode is uniquely attributed to one sample. This is a great tool to identify mistakes during index sequencing. Indeed, a phenomenon, known as index hopping, can lead to misattribution of some reads to the wrong sample. This is particularly frequent with the NovaSeq6000, and thus the use of Unique Dual-Indexing (UDI) is therefore highly recommended when using this sequencer.

The 24 UDI for tagmented libraries – Set I includes 24 primer pairs for unique dual-indexing allowing the multiplexing of up to 24 samples for sequencing on Illumina platforms. This set of indexes is designed to be used with  $\mu$ ChIPmentation for Histones (Cat. No. C01011011) or TAG Kit for ChIPmentation (Cat. No. C01011030) but it is also compatible with other tagmentation-based library preparation protocols, such as ATAC-Seq or CUT&Tag technologies.

The primer indexes for tagmented libraries contain 8-nucleotide long i7 and i5 indexes.

Table 1. Components supplied with the 24 UDI for tagmented libraries - Set I and their indexes sequences

Component	i7 Index Sequence	i5 Index Sequence	Cap color	Volume	Storage
Primer pair UDI 1	AACCGCGG	AGCGCTAG	transparent	12 $\mu$ l	-20°C
Primer pair UDI 2	GGTTATAA	GATATCGA	none	5 $\mu$ l	-20°C
Primer pair UDI 3	CCAAGTCC	CGCAGACG	none	5 $\mu$ l	-20°C
Primer pair UDI 4	TTGGACTT	TATGAGTA	none	5 $\mu$ l	-20°C
Primer pair UDI 5	CAGTGGAT	AGGTGCGT	none	5 $\mu$ l	-20°C
Primer pair UDI 6	TGACAAGC	GAACATAC	none	5 $\mu$ l	-20°C
Primer pair UDI 7	CTAGCTTG	ACATAGCG	none	5 $\mu$ l	-20°C

Component	i7 Index Sequence	i5 Index Sequence	Cap color	Volume	Storage
Primer pair UDI 8	TCGATCCA	GTGCGATA	none	5 µl	-20°C
Primer pair UDI 9	CCTGAACT	CCAACAGA	none	5 µl	-20°C
Primer pair UDI 10	TTCAGGTC	TTGGTGAG	none	5 µl	-20°C
Primer pair UDI 11	AGTAGAGA	CGCGGTTC	none	5 µl	-20°C
Primer pair UDI 12	GACGAGAG	TATAACCT	none	5 µl	-20°C
Primer pair UDI 13	AGACTTGG	AAGGATGA	none	5 µl	-20°C
Primer pair UDI 14	GAGTCCAA	GGAAGCAG	none	5 µl	-20°C
Primer pair UDI 15	CTTAAGCC	TCGTGACC	none	5 µl	-20°C
Primer pair UDI 16	TCCGGATT	CTACAGTT	none	5 µl	-20°C
Primer pair UDI 17	CTGTATTA	ATATTCAC	none	5 µl	-20°C
Primer pair UDI 18	TCACGCCG	GCGCCTGT	none	5 µl	-20°C
Primer pair UDI 19	ACTTACAT	ACTCTATG	none	5 µl	-20°C
Primer pair UDI 20	GTCCGTGC	GTCTCGCA	none	5 µl	-20°C
Primer pair UDI 21	AAGGTACC	AAGACGTC	none	5 µl	-20°C
Primer pair UDI 22	GGAACGTT	GGAGTACT	none	5 µl	-20°C
Primer pair UDI 23	AATTCTGC	ACCGGCCA	none	5 µl	-20°C
Primer pair UDI 24	GGCCTCAT	GTTAATTG	none	5 µl	-20°C

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

# 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	UDI 1 – UDI 10 UDI 2 – UDI 9 UDI 3 – UDI 12 UDI 4 – UDI 11 UDI 5 – UDI 14 UDI 6 – UDI 13 UDI 7 – UDI 16 UDI 8 – UDI 15
3 samples	UDI 1 – UDI 2 – UDI 9 UDI 3 – UDI 4 – UDI 11 UDI 5 – UDI 6 – UDI 13 UDI 7 – UDI 8 – UDI 15
4 samples	UDI 1 – UDI 2 – UDI 9 – UDI 10 UDI 3 – UDI 4 – UDI 11 – UDI 12 UDI 5 – UDI 6 – UDI 13 – UDI 14 UDI 7 – UDI 8 – UDI 15 – UDI 16

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.

# 24 UDI for tagmented libraries - Set II

Cat. No. C01011036

The 24 UDI (Unique dual indexes) for tagmented libraries – Set II provides combinations of barcodes where each barcode is uniquely attributed to one sample. This is a great tool to identify mistakes during index sequencing. Indeed, a phenomenon, known as index hopping, can lead to misattribution of some reads to the wrong sample. This is particularly frequent with the NovaSeq6000, and thus the use of Unique Dual-Indexing (UDI) is therefore highly recommended when using this sequencer.

The 24 UDI for tagmented libraries – Set II includes 24 primer pairs for unique dual-indexing and is designed to be combined with the 24 UDI for tagmented libraries – Set I, allowing multiplexing of up to 48 samples for sequencing on Illumina platforms. Both sets of indexes are designed to be used with ChIPmentation Kit for Histones (Cat. No. C01011009),  $\mu$ ChIPmentation for Histones (Cat. No. C01011011) or TAG Kit for ChIPmentation (Cat. No. C01011030) but it is also compatible with other tagmentation-based library preparation protocols, such as ATAC-Seq or CUT&Tag technologies. The primer indexes for tagmented libraries contain 8-nucleotide long i7 and i5 indexes.

**Table 1.** Components supplied with the 24 UDI for tagmented libraries – Set II and their indexes sequences

Component	i7 Index Sequence	i5 Index Sequence	Cap color	Volume	Storage
Primer pair UDI 25	ATCTTAGT	AACCGCGG	none	5 $\mu$ l	-20°C
Primer pair UDI 26	GCTCCGAC	GGTTATAA	none	5 $\mu$ l	-20°C
Primer pair UDI 27	ATACCAAG	CCAAGTCC	none	5 $\mu$ l	-20°C
Primer pair UDI 28	GCGTTGGA	TTGGACTT	none	5 $\mu$ l	-20°C
Primer pair UDI 29	CTTCACGG	CAGTGGAT	none	5 $\mu$ l	-20°C
Primer pair UDI 30	TCCTGTAA	TGACAAGC	none	5 $\mu$ l	-20°C
Primer pair UDI 31	AGAATGCC	CTAGCTTG	none	5 $\mu$ l	-20°C



Component	i7 Index Sequence	i5 Index Sequence	Cap color	Volume	Storage
Primer pair UDI 32	GAGGCATT	TCGATCCA	none	5 µl	-20°C
Primer pair UDI 33	CCTCGGTA	CCTGAACT	none	5 µl	-20°C
Primer pair UDI 34	TTCTAACG	TTCAGGTC	none	5 µl	-20°C
Primer pair UDI 35	ATGAGGCT	AGTAGAGA	none	5 µl	-20°C
Primer pair UDI 36	GCAGAATC	GACGAGAG	none	5 µl	-20°C
Primer pair UDI 37	CACTACGA	AGACTTGG	none	5 µl	-20°C
Primer pair UDI 38	TGTCGTAG	GAGTCCAA	none	5 µl	-20°C
Primer pair UDI 39	ACCACTTA	CTTAAGCC	none	5 µl	-20°C
Primer pair UDI 40	GTTGTCCG	TCCGGATT	none	5 µl	-20°C
Primer pair UDI 41	ATCCATAT	CTGTATTA	none	5 µl	-20°C
Primer pair UDI 42	GCTTGCGC	TCACGCCG	none	5 µl	-20°C
Primer pair UDI 43	AGTATCTT	ACTTACAT	none	5 µl	-20°C
Primer pair UDI 44	GACGCTCC	GTCCGTGC	none	5 µl	-20°C
Primer pair UDI 45	CATGCCAT	AAGGTACC	none	5 µl	-20°C
Primer pair UDI 46	TGCATTGC	GGAACGTT	none	5 µl	-20°C
Primer pair UDI 47	ATTGGAAC	AATTCTGC	none	5 µl	-20°C
Primer pair UDI 48	GCCAAGGT	GGCCTCAT	none	5 µl	-20°C

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

# 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	UDI 17 – UDI 26 UDI 18 – UDI 25 UDI 19 – UDI 28 UDI 20 – UDI 27 UDI 21 – UDI 30 UDI 22 – UDI 29 UDI 23 – UDI 32 UDI 24 – UDI 31 UDI 33 – UDI 42 UDI 34 – UDI 41 UDI 35 – UDI 44 UDI 36 – UDI 43 UDI 37 – UDI 46 UDI 38 – UDI 45 UDI 39 – UDI 48 UDI 40 – UDI 47
3 samples	UDI 17 – UDI 18 – UDI 25 UDI 19 – UDI 20 – UDI 27 UDI 21 – UDI 22 – UDI 29 UDI 23 – UDI 24 – UDI 31 UDI 33 – UDI 34 – UDI 41 UDI 35 – UDI 36 – UDI 43 UDI 37 – UDI 38 – UDI 45 UDI 39 – UDI 40 – UDI 47
4 samples	UDI 17 – UDI 18 – UDI 25 – UDI 26 UDI 19 – UDI 20 – UDI 27 – UDI 28 UDI 21 – UDI 22 – UDI 29 – UDI 29 UDI 23 – UDI 24 – UDI 31 – UDI 32 UDI 33 – UDI 34 – UDI 41 – UDI 42 UDI 35 – UDI 36 – UDI 43 – UDI 44 UDI 37 – UDI 38 – UDI 45 – UDI 46 UDI 39 – UDI 40 – UDI 47 – UDI 48

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.
ChIPmentation Kit for Histones (24 rxns)	C01011009
$\mu$ ChIPmentation Kit for Histones (24 rxns)	C01011011
$\mu$ ChIPmentation Kit for Histones (8 rxns)	C01011012
TAG Kit for ChIPmentation (24 rxns)	C01011030
Tagmentase (loaded)	C01070012
Tagmentase	C01070010
Tagmentase Dilution Buffer	C01070011
pA-Tn5 Transposase - loaded	C01070001
pA-Tn5 Transposase - unloaded	C01070002

## **FOR RESEARCH USE ONLY.**

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

© 2020 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)