

# Recommended Alternative Microarray Consumables

Pub. No. MAN0019853 Rev. H00

## Contents

- “Axiom™ assays manual workflows” on page 1.
- “Axiom™ assays automated workflows—NIMBUS® Target Preparation Instrument” on page 2.
- “Axiom™ assays automated workflows—Beckman Coulter™ Biomek™ liquid handlers” on page 3.
- “Axiom™ PharmacoFocus™ assays” on page 3.
- “Axiom™ 2.0 PLUS assays” on page 4.
- “CarrierScan™ assays” on page 5.
- “CytoScan™ assays” on page 6.
- “Expression assays” on page 7.
- “PharmacoScan™ assays” on page 7.
- “Other kits” on page 9.

## Overview

In the event of plate shortages from suppliers, alternative consumables that have been tested in the laboratories of Thermo Fisher Scientific can be directly and safely substituted. The following tables list the tested and recommended alternative consumables for the assays described.

**IMPORTANT!** For any consumable item shortage not listed in this document, contact your local FAS to help manage the issue. Use of other untested consumables can result in assay failure and may violate the array and reagent replacement policy.

Unless otherwise indicated, all materials are available through [thermofisher.com](https://www.thermofisher.com).

## Axiom™ assays manual workflows

| Consumable                                                        |                                   |                   | Recommended replacement                                                         |                                   |                                                    |
|-------------------------------------------------------------------|-----------------------------------|-------------------|---------------------------------------------------------------------------------|-----------------------------------|----------------------------------------------------|
| Name                                                              | Component Part No. <sup>[1]</sup> | Source            | Name                                                                            | Component Part No. <sup>[1]</sup> | Source                                             |
| <b>Axiom™ 2.0 Assay 24-Array Format Manual Workflow</b>           |                                   |                   |                                                                                 |                                   |                                                    |
| Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (blue)   | N/A                               | Bio-Rad™, HSP9631 | Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (white) <sup>[2]</sup> | N/A                               | Bio-Rad™, HSP9601                                  |
| Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (green) | N/A                               | Bio-Rad™, HSS9641 | MicroAmp™ EnduraPlate™ Optical 96-Well Clear Reaction Plates with Barcode       | N/A                               | <a href="#">4483354</a> or <a href="#">4483352</a> |
|                                                                   |                                   |                   | Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (clear)               | N/A                               | Bio-Rad™, HSS9601                                  |

| Consumable                                                         |                                   |                                                | Recommended replacement                                                         |                                   |                                                    |
|--------------------------------------------------------------------|-----------------------------------|------------------------------------------------|---------------------------------------------------------------------------------|-----------------------------------|----------------------------------------------------|
| Name                                                               | Component Part No. <sup>[1]</sup> | Source                                         | Name                                                                            | Component Part No. <sup>[1]</sup> | Source                                             |
| Axiom™ 2.0 Assay 96-Array Format Manual Workflow                   |                                   |                                                |                                                                                 |                                   |                                                    |
| Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (blue)    | N/A                               | Bio-Rad™, HSP9631                              | Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (white) <sup>[2]</sup> | N/A                               | Bio-Rad™, HSP9601                                  |
| Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (green)  | N/A                               | Bio-Rad™, HSS9641                              | MicroAmp™ EnduraPlate™ Optical 96-Well Clear Reaction Plates with Barcode       | N/A                               | <a href="#">4483354</a> or <a href="#">4483352</a> |
|                                                                    |                                   |                                                | Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (clear)               | N/A                               | Bio-Rad™, HSS9601                                  |
| Axiom™ 2.0 Assay Mini 96-Array Format Manual Workflow              |                                   |                                                |                                                                                 |                                   |                                                    |
| Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (blue)    | 203015                            | Bio-Rad™, HSP9631                              | Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (white) <sup>[2]</sup> | 203208                            | Bio-Rad™, HSP9601                                  |
| Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (green)  | 203009                            | Bio-Rad™, HSS9641                              | MicroAmp™ EnduraPlate™ Optical 96-Well Clear Reaction Plates with Barcode       | 203244                            | <a href="#">4483354</a> or <a href="#">4483352</a> |
| Eppendorf™ DeepWell™ Plate 96 ( <a href="#">951033481</a> )        | 203079                            | Fisher Scientific™, <a href="#">13-864-302</a> | Thermo Scientific™ Abgene™ 96 Well 2.2 mL Polypropylene Deepwell Storage Plate  | 202920                            | Fisher Scientific™, <a href="#">AB0932</a>         |
| Axiom™ 2.0 Assay Mini 96 Manual Target Preparation Consumables Kit |                                   | <a href="#">902986</a>                         | Axiom™ 2.0 Assay Mini 96 Manual Target Preparation Consumables Kit v2           |                                   | <a href="#">952431</a>                             |

<sup>[1]</sup> As a part of a larger Applied Biosystems™ consumables kit, the component part number is used for identification purposes only, and is not an orderable catalog number.

<sup>[2]</sup> The Hard-Shell™ 96-Well PCR Plate is available in several colors, any of which are acceptable.

## Axiom™ assays automated workflows—NIMBUS® Target Preparation Instrument

| Consumable                                                                                        |                                   |                                                | Recommended replacement                                                            |                                   |                                                    |
|---------------------------------------------------------------------------------------------------|-----------------------------------|------------------------------------------------|------------------------------------------------------------------------------------|-----------------------------------|----------------------------------------------------|
| Name                                                                                              | Component Part No. <sup>[1]</sup> | Source                                         | Name                                                                               | Component Part No. <sup>[1]</sup> | Source                                             |
| Axiom™ 2.0 Assay 96-Array Format Automated Workflow on the Applied Biosystems™ NIMBUS® Instrument |                                   |                                                |                                                                                    |                                   |                                                    |
| Round deepwell plate<br>Axygen™ 96-well Round Deepwell Plate (sterile, PDW20CS)                   | 203028                            | Fisher Scientific™, <a href="#">14-222-354</a> | Round deepwell plate<br>Axygen™ 96-well Round Deepwell Plate (non-sterile, PDW20C) | 203367                            | Fisher Scientific™, <a href="#">14-222-353</a>     |
| Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (green)                                 | 203009                            | Bio-Rad™, HSS9641                              | MicroAmp™ EnduraPlate™ Optical 96-Well Clear Reaction Plates with Barcode          | 203244                            | <a href="#">4483354</a> or <a href="#">4483352</a> |
|                                                                                                   |                                   |                                                | Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (clear)                  | N/A                               | Bio-Rad™, HSS9601                                  |
| Axiom™ 96 Consumables Kit for Applied Biosystems™ NIMBUS® 2.0                                     |                                   | <a href="#">902907</a>                         | Axiom™ NIMBUS® 2.0 Consumables Kit v2                                              |                                   | <a href="#">952437</a>                             |

<sup>[1]</sup> As a part of a larger Applied Biosystems™ consumables kit, the component part number is used for identification purposes only, and is not an orderable catalog number.

## Axiom™ assays automated workflows—Beckman Coulter™ Biomek™ liquid handlers

| Consumable                                                                                                                                                                                  |                                   |                                                | Recommended replacement                                                         |                                   |                                            |
|---------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|-----------------------------------|------------------------------------------------|---------------------------------------------------------------------------------|-----------------------------------|--------------------------------------------|
| Name                                                                                                                                                                                        | Component Part No. <sup>[1]</sup> | Source                                         | Name                                                                            | Component Part No. <sup>[1]</sup> | Source                                     |
| <b>Axiom™ 2.0 Assay 96-Array Format Automated Workflow on the Biomek™ FX<sup>P</sup> Target Prep Express Automated Workstation</b>                                                          |                                   |                                                |                                                                                 |                                   |                                            |
| Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (blue)                                                                                                                             | 203015                            | Bio-Rad™, HSP9631                              | Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (white) <sup>[2]</sup> | 203208                            | Bio-Rad™, HSP9601                          |
| Axiom™ 96-format Consumables Kit for QC (Windows™ 7)                                                                                                                                        | Greiner, EK-25801                 | 902801                                         | Axiom™ Consumables Kit for QC                                                   | Corning, 3635                     | <a href="#">902909</a>                     |
| <b>Axiom™ 2.0 Assay 384HT Array Format Automated Workflow on the Biomek™ i7</b>                                                                                                             |                                   |                                                |                                                                                 |                                   |                                            |
| Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (blue)                                                                                                                             | 203015                            | Bio-Rad™, HSP9631                              | Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (white) <sup>[2]</sup> | 203208                            | Bio-Rad™, HSP9601                          |
| Axygen™ 384-well Clear 240 µL Polypropylene Deep Well Microplate (non-barcoded)                                                                                                             | 203103                            | Fisher Scientific™, <a href="#">14-222-229</a> | Abgene™ 384-Well 250 µL Polypropylene Storage Plate (non-barcoded)              | 203251                            | Fisher Scientific™, <a href="#">AB1178</a> |
| Axygen™ 384-well Clear 240 µL Polypropylene Deep Well Microplate (barcoded)                                                                                                                 | 203017                            | Fisher Scientific™, <a href="#">14-222-229</a> | Abgene™ 384-Well 250 µL Polypropylene Storage Plate (barcoded)                  | 203294                            | Fisher Scientific™, <a href="#">AB1178</a> |
| <b>Axiom™ 2.0 Assay 384HT Array Format Automated Workflow on the Biomek™ FX<sup>P</sup> Target Prep Express Automated Workstation for Windows™ XP and Windows™ 7 versions of the method</b> |                                   |                                                |                                                                                 |                                   |                                            |
| Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (blue)                                                                                                                             | 203015                            | Bio-Rad™, HSP9631                              | Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (white) <sup>[2]</sup> | 203208                            | Bio-Rad™, HSP9601                          |
| Axiom™ 384HT Consumables Kit for Biomek™ FX <sup>P</sup>                                                                                                                                    |                                   | <a href="#">902288</a>                         | Axiom™ 384HT Consumables Kit for Biomek™ FX <sup>P</sup> v2                     |                                   | <a href="#">952455</a> <sup>[3]</sup>      |

<sup>[1]</sup> As a part of a larger Applied Biosystems™ consumables kit, the component part number is used for identification purposes only, and is not an orderable catalog number.

<sup>[2]</sup> The Hard-Shell™ 96-Well PCR Plate is available in several colors, any of which are acceptable.

<sup>[3]</sup> Axiom™ 384HT Consumables Kit for Biomek™ FX<sup>P</sup> ([902288](#)), contains 15 x 203017 (barcoded). Axiom™ 384HT Consumables Kit for Biomek™ FX<sup>P</sup> v2 ([952455](#)) contains 10 x 203017 (barcoded) and 5 x 203251 (non-barcoded) for a total of 15 x 384 deepwell plates.

## Axiom™ PharmacoFocus™ assays

| Consumable                                                              |                                   |                   | Recommended replacement                                                         |                                   |                                                    |
|-------------------------------------------------------------------------|-----------------------------------|-------------------|---------------------------------------------------------------------------------|-----------------------------------|----------------------------------------------------|
| Name                                                                    | Component Part No. <sup>[1]</sup> | Source            | Name                                                                            | Component Part No. <sup>[1]</sup> | Source                                             |
| <b>Axiom™ PharmacoFocus™ Assay Mini 96-Array Format Manual Workflow</b> |                                   |                   |                                                                                 |                                   |                                                    |
| Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (blue)         | N/A                               | Bio-Rad™, HSP9631 | Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (white) <sup>[2]</sup> | N/A                               | Bio-Rad™, HSP9601                                  |
| Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (green)       | N/A                               | Bio-Rad™, HSS9641 | MicroAmp™ EnduraPlate™ Optical 96-Well Clear Reaction Plates with Barcode       | N/A                               | <a href="#">4483354</a> or <a href="#">4483352</a> |
|                                                                         |                                   |                   | Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (clear)               | N/A                               | Bio-Rad™, HSS9601                                  |

| Consumable                                                                                                               |                                   |                                                | Recommended replacement                                                                      |                                   |                                                    |
|--------------------------------------------------------------------------------------------------------------------------|-----------------------------------|------------------------------------------------|----------------------------------------------------------------------------------------------|-----------------------------------|----------------------------------------------------|
| Name                                                                                                                     | Component Part No. <sup>[1]</sup> | Source                                         | Name                                                                                         | Component Part No. <sup>[1]</sup> | Source                                             |
| <b>Axiom™ PharmacoFocus™ Assay Mini 96-Array Format Automated Workflow on the Applied Biosystems™ NIMBUS® Instrument</b> |                                   |                                                |                                                                                              |                                   |                                                    |
| Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (blue) for mPCR use                                             | 203015                            | Bio-Rad™, HSP9631                              | Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (white) for mPCR use <sup>[2]</sup> | 203023                            | Bio-Rad™, HSP9601                                  |
| Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (green) for mPCR use                                           | 203009                            | Bio-Rad™, HSS9641                              | MicroAmp™ EnduraPlate™ Optical 96-Well Clear Reaction Plates with Barcode for mPCR use       | 203244                            | <a href="#">4483354</a> or <a href="#">4483352</a> |
|                                                                                                                          |                                   |                                                | Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (clear) for mPCR use               | N/A                               | Bio-Rad™, HSS9601                                  |
| Round deepwell plate<br>Axygen™ 96-well Round Deepwell Plate (sterile, PDW20CS)                                          | 203028                            | Fisher Scientific™, <a href="#">14-222-354</a> | Round deepwell plate<br>Axygen™ 96-well Round Deepwell Plate (non-sterile, PDW20C)           | 203367                            | Fisher Scientific™, <a href="#">14-222-353</a>     |
| Axiom™ 96 Consumables Kit for Applied Biosystems™ NIMBUS® 2.0                                                            |                                   | <a href="#">902907</a>                         | Axiom™ NIMBUS® 2.0 Consumables Kit v2                                                        |                                   | <a href="#">952437</a>                             |

<sup>[1]</sup> As a part of a larger Applied Biosystems™ consumables kit, the component part number is used for identification purposes only, and is not an orderable catalog number.

## Axiom™ 2.0 PLUS assays

| Consumable                                                                                                    |                                   |                                                | Recommended replacement                                                                      |                                   |                                                    |
|---------------------------------------------------------------------------------------------------------------|-----------------------------------|------------------------------------------------|----------------------------------------------------------------------------------------------|-----------------------------------|----------------------------------------------------|
| Name                                                                                                          | Component Part No. <sup>[1]</sup> | Source                                         | Name                                                                                         | Component Part No. <sup>[1]</sup> | Source                                             |
| <b>Axiom™ 2.0 Plus Assay 96-Array Format Automated Workflow on the Applied Biosystems™ NIMBUS® Instrument</b> |                                   |                                                |                                                                                              |                                   |                                                    |
| Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (blue) for mPCR use                                  | 203015                            | Bio-Rad™, HSP9631                              | Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (white) for mPCR use <sup>[2]</sup> | 203023                            | Bio-Rad™, HSP9601                                  |
| Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (green) for mPCR use                                | 203009                            | Bio-Rad™, HSS9641                              | MicroAmp™ EnduraPlate™ Optical 96-Well Clear Reaction Plates with Barcode for mPCR use       | 203244                            | <a href="#">4483354</a> or <a href="#">4483352</a> |
|                                                                                                               |                                   |                                                | Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (clear) for mPCR use               | N/A                               | Bio-Rad™, HSS9601                                  |
| Round deepwell plate<br>Axygen™ 96-well Round Deepwell Plate (sterile, PDW20CS)                               | 203028                            | Fisher Scientific™, <a href="#">14-222-354</a> | Round deepwell plate<br>Axygen™ 96-well Round Deepwell Plate (non-sterile, PDW20C)           | 203367                            | Fisher Scientific™, <a href="#">14-222-353</a>     |
| Axiom™ 96 Consumables Kit for Applied Biosystems™ NIMBUS® 2.0                                                 |                                   | <a href="#">902907</a>                         | Axiom™ NIMBUS® 2.0 Consumables Kit v2                                                        |                                   | <a href="#">952437</a>                             |

<sup>[1]</sup> As a part of a larger Applied Biosystems™ consumables kit, the component part number is used for identification purposes only, and is not an orderable catalog number.

<sup>[2]</sup> The Hard-Shell™ 96-Well PCR Plate is available in several colors, any of which are acceptable.

## CarrierScan™ assays

| Consumable                                                                                                 |                                   |                                                | Recommended replacement                                                                      |                                   |                                                    |
|------------------------------------------------------------------------------------------------------------|-----------------------------------|------------------------------------------------|----------------------------------------------------------------------------------------------|-----------------------------------|----------------------------------------------------|
| Name                                                                                                       | Component Part No. <sup>[1]</sup> | Source                                         | Name                                                                                         | Component Part No. <sup>[1]</sup> | Source                                             |
| <b>CarrierScan™ Assay 96-Array Format Manual Workflow</b>                                                  |                                   |                                                |                                                                                              |                                   |                                                    |
| Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (blue)                                            | N/A                               | Bio-Rad™, HSP9631                              | Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (white) <sup>[2]</sup>              | N/A                               | Bio-Rad™, HSP9601                                  |
| Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (green)                                          | N/A                               | Bio-Rad™, HSS9641                              | MicroAmp™ EnduraPlate™ Optical 96-Well Clear Reaction Plates with Barcode                    | N/A                               | <a href="#">4483354</a> or <a href="#">4483352</a> |
|                                                                                                            |                                   |                                                | Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (clear)                            | N/A                               | Bio-Rad™, HSS9601                                  |
| Eppendorf™ DeepWell™ Plate 96 ( <a href="#">951033481</a> )                                                | N/A                               | Fisher Scientific™, <a href="#">13-864-302</a> | Thermo Scientific™ Abgene™ 96 Well 2.2 mL Polypropylene Deepwell Storage Plate               | N/A                               | Fisher Scientific™, <a href="#">AB0932</a>         |
| Axiom™ 96-format Consumables Kit for QC (Windows™ 7)                                                       | Greiner, EK-25801                 | 902801                                         | Axiom™ Consumables Kit for QC                                                                | Corning, 3635                     | <a href="#">902909</a>                             |
| <b>CarrierScan™ Assay 96-Array Format Automated Workflow on the Applied Biosystems™ NIMBUS® Instrument</b> |                                   |                                                |                                                                                              |                                   |                                                    |
| Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (blue) for mPCR use                               | N/A                               | Bio-Rad™, HSP9631                              | Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (white) for mPCR use <sup>[2]</sup> | N/A                               | Bio-Rad™, HSP9601                                  |
| Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (green) for mPCR use                             | N/A                               | Bio-Rad™, HSS9641                              | MicroAmp™ EnduraPlate™ Optical 96-Well Clear Reaction Plates with Barcode for mPCR use       | N/A                               | <a href="#">4483354</a> or <a href="#">4483352</a> |
|                                                                                                            |                                   |                                                | Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (clear) for mPCR use               | N/A                               | Bio-Rad™, HSS9601                                  |
| Round deepwell plate<br>Axygen™ 96-well Round Deepwell Plate (sterile, PDW20CS)                            | 203028                            | Fisher Scientific™, <a href="#">14-222-354</a> | Round deepwell plate<br>Axygen™ 96-well Round Deepwell Plate (non-sterile, PDW20C)           | 203367                            | Fisher Scientific™, <a href="#">14-222-353</a>     |
| HT Target Prep 96 Consumables Kit for Applied Biosystems™ NIMBUS® 2.0                                      |                                   | <a href="#">951996</a>                         | HT Target Prep 96 Consumables Kit for Applied Biosystems™ NIMBUS® 2.0, v2                    |                                   | <a href="#">952438</a>                             |

| Consumable                                                                                                                           |                                   |                                                | Recommended replacement                                                                      |                                   |                                                    |
|--------------------------------------------------------------------------------------------------------------------------------------|-----------------------------------|------------------------------------------------|----------------------------------------------------------------------------------------------|-----------------------------------|----------------------------------------------------|
| Name                                                                                                                                 | Component Part No. <sup>[1]</sup> | Source                                         | Name                                                                                         | Component Part No. <sup>[1]</sup> | Source                                             |
| <b>CarrierScan™ Assay 96-Array Format Automated Workflow on the Biomek™ FX<sup>P</sup> Target Prep Express Automated Workstation</b> |                                   |                                                |                                                                                              |                                   |                                                    |
| Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (blue) for mPCR use                                                         | N/A                               | Bio-Rad™, HSP9631                              | Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (white) for mPCR use <sup>[2]</sup> | N/A                               | Bio-Rad™, HSP9601                                  |
| Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (green) for mPCR use                                                       | N/A                               | Bio-Rad™, HSS9641                              | MicroAmp™ EnduraPlate™ Optical 96-Well Clear Reaction Plates with Barcode for mPCR use       | N/A                               | <a href="#">4483354</a> or <a href="#">4483352</a> |
|                                                                                                                                      |                                   |                                                | Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (clear) for mPCR use               | N/A                               | Bio-Rad™, HSS9601                                  |
| Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (blue)                                                                      | 203015                            | Bio-Rad™, HSP9631                              | Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (white) <sup>[2]</sup>              | 203208                            | Bio-Rad™, HSP9601                                  |
| Eppendorf™ DeepWell™ Plate 96 ( <a href="#">951033481</a> )                                                                          | 203079                            | Fisher Scientific™, <a href="#">13-864-302</a> | Thermo Scientific™ Abgene™ 96 Well 2.2 mL Polypropylene Deepwell Storage Plate               | 202920                            | Fisher Scientific™, <a href="#">AB0932</a>         |
| Axiom™ 96-format Consumables Kit for Biomek™ FX <sup>P</sup> (Windows™ 7)                                                            |                                   | 902800                                         | Axiom™ 96-Format Consumables Kit for Biomek™ FX <sup>P</sup> Windows™ 7, v2                  |                                   | 952430                                             |
| Axiom™ 96-format Consumables Kit for QC (Windows™ 7)                                                                                 | Greiner, EK-25801                 | 902801                                         | Axiom™ Consumables Kit for QC                                                                | Corning, 3635                     | <a href="#">902909</a>                             |

<sup>[1]</sup> As a part of a larger Applied Biosystems™ consumables kit, the component part number is used for identification purposes only, and is not an orderable catalog number.

<sup>[2]</sup> The Hard-Shell™ 96-Well PCR Plate is available in several colors, any of which are acceptable.

## CytoScan™ assays

| Consumable                                                        |                                   |                                                | Recommended replacement                                                         |                                   |                                                    |
|-------------------------------------------------------------------|-----------------------------------|------------------------------------------------|---------------------------------------------------------------------------------|-----------------------------------|----------------------------------------------------|
| Name                                                              | Component Part No. <sup>[1]</sup> | Source                                         | Name                                                                            | Component Part No. <sup>[1]</sup> | Source                                             |
| <b>CytoScan™ HT-CMA Assay 96-Array Format Manual Workflow</b>     |                                   |                                                |                                                                                 |                                   |                                                    |
| Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (blue)   | N/A                               | Bio-Rad™, HSP9631                              | Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (white) <sup>[2]</sup> | N/A                               | Bio-Rad™, HSP9601                                  |
| Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (green) | N/A                               | Bio-Rad™, HSS9641                              | MicroAmp™ EnduraPlate™ Optical 96-Well Clear Reaction Plates with Barcode       | N/A                               | <a href="#">4483354</a> or <a href="#">4483352</a> |
|                                                                   |                                   |                                                | Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (clear)               | N/A                               | Bio-Rad™, HSS9601                                  |
| Eppendorf™ DeepWell™ Plate 96 ( <a href="#">951033481</a> )       | N/A                               | Fisher Scientific™, <a href="#">13-864-302</a> | Thermo Scientific™ Abgene™ 96 Well 2.2 mL Polypropylene Deepwell Storage Plate  | N/A                               | Fisher Scientific™, <a href="#">AB0932</a>         |

| Consumable                                                                                              |                                   |                                                | Recommended replacement                                                                                                                                                                                   |                                   |                                                                                                                                                        |
|---------------------------------------------------------------------------------------------------------|-----------------------------------|------------------------------------------------|-----------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|-----------------------------------|--------------------------------------------------------------------------------------------------------------------------------------------------------|
| Name                                                                                                    | Component Part No. <sup>[1]</sup> | Source                                         | Name                                                                                                                                                                                                      | Component Part No. <sup>[1]</sup> | Source                                                                                                                                                 |
| CytoScan™ HT-CMA Assay 96-Array Format Automated Workflow on the Applied Biosystems™ NIMBUS® Instrument |                                   |                                                |                                                                                                                                                                                                           |                                   |                                                                                                                                                        |
| Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (green)                                       | 203009                            | Bio-Rad™, HSS9641                              | Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (clear)                                                                                                                                         | N/A                               | Bio-Rad™, HSS9601                                                                                                                                      |
| Round deepwell plate<br>Axygen™ 96-well Round Deepwell Plate (sterile, PDW20CS)                         | 203028                            | Fisher Scientific™, <a href="#">14-222-354</a> | Round deepwell plate <ul style="list-style-type: none"><li>Axygen™ 96-well Round Deepwell Plate (non-sterile, PDW20C)</li><li>PerkinElmer™ Deepwell StorPlate-96 U Round, 2.0 mL (6008490 only)</li></ul> | 203367                            | <ul style="list-style-type: none"><li>Fisher Scientific™, <a href="#">14-222-353</a></li><li>Fisher Scientific™, <a href="#">50-905-1741</a></li></ul> |
| HT Target Prep 96 Consumables Kit for Applied Biosystems™ NIMBUS® 2.0                                   |                                   | <a href="#">951996</a>                         | HT Target Prep 96 Consumables Kit for Applied Biosystems™ NIMBUS® 2.0, v2                                                                                                                                 |                                   | <a href="#">952438</a>                                                                                                                                 |
| CytoScan™ Assay Automated Workflow on the Applied Biosystems™ NIMBUS® Instrument (25 or 49 samples)     |                                   |                                                |                                                                                                                                                                                                           |                                   |                                                                                                                                                        |
| Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (green)                                       | 203009                            | Bio-Rad™, HSS9641                              | MicroAmp™ EnduraPlate™ Optical 96-Well Clear Reaction Plates with Barcode                                                                                                                                 | 203244                            | <a href="#">4483354</a> or <a href="#">4483352</a>                                                                                                     |
| CytoScan™ Consumables Kit for NIMBUS®                                                                   |                                   | <a href="#">902168</a>                         | CytoScan™ Consumables Kit for NIMBUS® v2                                                                                                                                                                  |                                   | <a href="#">952449</a>                                                                                                                                 |

<sup>[1]</sup> As a part of a larger Applied Biosystems™ consumables kit, the component part number is used for identification purposes only, and is not an orderable catalog number.

<sup>[2]</sup> The Hard-Shell™ 96-Well PCR Plate is available in several colors, any of which are acceptable.

## Expression assays

| Consumable                                                      |                                   |                   | Recommended replacement                                                         |                                   |                   |
|-----------------------------------------------------------------|-----------------------------------|-------------------|---------------------------------------------------------------------------------|-----------------------------------|-------------------|
| Name                                                            | Component Part No. <sup>[1]</sup> | Source            | Name                                                                            | Component Part No. <sup>[1]</sup> | Source            |
| <b>Expression microarray assays (all)</b>                       |                                   |                   |                                                                                 |                                   |                   |
| Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (blue) | 202916                            | Bio-Rad™, HSP9631 | Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (white) <sup>[2]</sup> | 203208                            | Bio-Rad™, HSP9601 |
| Plate, UV Star                                                  | 202166                            | Greiner EK-25801  | Plate, UV Transparent Flat 96 wells                                             | 202919                            | Corning 3625      |

<sup>[1]</sup> As a part of a larger Applied Biosystems™ consumables kit, the component part number is used for identification purposes only, and is not an orderable catalog number.

<sup>[2]</sup> The Hard-Shell™ 96-Well PCR Plate is available in several colors, any of which are acceptable.

## PharmacoScan™ assays

| Consumable                                                        |                                   |                   | Recommended replacement                                                         |                                   |                                                    |
|-------------------------------------------------------------------|-----------------------------------|-------------------|---------------------------------------------------------------------------------|-----------------------------------|----------------------------------------------------|
| Name                                                              | Component Part No. <sup>[1]</sup> | Source            | Name                                                                            | Component Part No. <sup>[1]</sup> | Source                                             |
| <b>PharmacoScan™ Assay 24-Array Format Manual Workflow</b>        |                                   |                   |                                                                                 |                                   |                                                    |
| Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (blue)   | N/A                               | Bio-Rad™, HSP9631 | Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (white) <sup>[2]</sup> | N/A                               | Bio-Rad™, HSP9601                                  |
| Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (green) | N/A                               | Bio-Rad™, HSS9641 | MicroAmp™ EnduraPlate™ Optical 96-Well Clear Reaction Plates with Barcode       | N/A                               | <a href="#">4483354</a> or <a href="#">4483352</a> |
|                                                                   |                                   |                   | Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (clear)               | N/A                               | Bio-Rad™, HSS9601                                  |

| Consumable                                                                                                                      |                                   |                                | Recommended replacement                                                                                                         |                                   |                                                                                |
|---------------------------------------------------------------------------------------------------------------------------------|-----------------------------------|--------------------------------|---------------------------------------------------------------------------------------------------------------------------------|-----------------------------------|--------------------------------------------------------------------------------|
| Name                                                                                                                            | Component Part No. <sup>[1]</sup> | Source                         | Name                                                                                                                            | Component Part No. <sup>[1]</sup> | Source                                                                         |
| PharmacoScan™ Assay 96-Array Format Manual Workflow                                                                             |                                   |                                |                                                                                                                                 |                                   |                                                                                |
| Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (blue)                                                                 | N/A                               | Bio-Rad™, HSP9631              | Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (white) <sup>[2]</sup>                                                 | N/A                               | Bio-Rad™, HSP9601                                                              |
| Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (green)                                                               | N/A                               | Bio-Rad™, HSS9641              | MicroAmp™ EnduraPlate™ Optical 96-Well Clear Reaction Plates with Barcode                                                       | N/A                               | 4483354 or 4483352                                                             |
|                                                                                                                                 |                                   |                                | Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (clear)                                                               | N/A                               | Bio-Rad™, HSS9601                                                              |
| PharmacoScan™ Assay 96-Array Format Automated Workflow on the Biomek™ FX <sup>P</sup> Target Prep Express Automated Workstation |                                   |                                |                                                                                                                                 |                                   |                                                                                |
| Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (blue) for mPCR use                                                    | N/A                               | Bio-Rad™, HSP9631              | Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (white) for mPCR use <sup>[2]</sup>                                    | N/A                               | Bio-Rad™, HSP9601                                                              |
| Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (green) for mPCR use                                                  | N/A                               | Bio-Rad™, HSS9641              | MicroAmp™ EnduraPlate™ Optical 96-Well Clear Reaction Plates with Barcode for mPCR use                                          | N/A                               | 4483354 or 4483352                                                             |
|                                                                                                                                 |                                   |                                | Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (clear) for mPCR use                                                  | N/A                               | Bio-Rad™, HSS9601                                                              |
| Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (blue)                                                                 | 203015                            | Bio-Rad™, HSP9631              | Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (white) <sup>[2]</sup>                                                 | 203208                            | Bio-Rad™, HSP9601                                                              |
| Eppendorf™ DeepWell™ Plate 96 (951033481)                                                                                       | 203079                            | Fisher Scientific™, 13-864-302 | Thermo Scientific™ Abgene™ 96 Well 2.2 mL Polypropylene Deepwell Storage Plate                                                  | 202920                            | Fisher Scientific™, AB0932                                                     |
| Axiom™ 96-format Consumables Kit for Biomek™ FX <sup>P</sup> (Windows™ 7)                                                       |                                   | 902800                         | Axiom™ 96-Format Consumables Kit for Biomek™ FX <sup>P</sup> Windows™ 7, v2                                                     |                                   | 952430                                                                         |
| PharmacoScan™ Assay 96-Array Format Automated Workflow on the Applied Biosystems™ NIMBUS® Instrument                            |                                   |                                |                                                                                                                                 |                                   |                                                                                |
| Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (blue) for mPCR use                                                    | 203015                            | Bio-Rad™, HSP9631              | Hard-Shell™ 96-Well PCR Plate, low profile, full skirted (white) for mPCR use <sup>[2]</sup>                                    | 203023                            | Bio-Rad™, HSP9601                                                              |
| Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (green) for mPCR use                                                  | 203009                            | Bio-Rad™, HSS9641              | MicroAmp™ EnduraPlate™ Optical 96-Well Clear Reaction Plates with Barcode for mPCR use                                          | 203244                            | 4483354 or 4483352                                                             |
|                                                                                                                                 |                                   |                                | Hard-Shell™ 96-Well PCR Plate, high profile, semi skirted (clear) for mPCR use                                                  | N/A                               | Bio-Rad™, HSS9601                                                              |
| Round deepwell plate<br>Axygen™ 96-well Round Deepwell Plate (sterile, PDW20CS)                                                 | 203028                            | Fisher Scientific™, 14-222-354 | Round deepwell plate <ul style="list-style-type: none"><li>Axygen™ 96-well Round Deepwell Plate (non-sterile, PDW20C)</li></ul> | 203367                            | <ul style="list-style-type: none"><li>Fisher Scientific™, 14-222-353</li></ul> |
| Axiom™ 96 Consumables Kit for Applied Biosystems™ NIMBUS® 2.0                                                                   |                                   | 902907                         | Axiom™ NIMBUS® 2.0 Consumables Kit v2                                                                                           |                                   | 952437                                                                         |

<sup>[1]</sup> As a part of a larger Applied Biosystems™ consumables kit, the component part number is used for identification purposes only, and is not an orderable catalog number.

<sup>[2]</sup> The Hard-Shell™ 96-Well PCR Plate is available in several colors, any of which are acceptable.

Other kits

| Consumable                                                 |        | Recommended replacement                                                   |        |
|------------------------------------------------------------|--------|---------------------------------------------------------------------------|--------|
| Name                                                       | Source | Name                                                                      | Source |
| Axiom™ Consumables Kit for Applied Biosystems™ NIMBUS® 1.0 | 902366 | Axiom™ Consumables Kit for Applied Biosystems™ NIMBUS® Instrument 1.0, v2 | 952436 |

Limited product warranty

Life Technologies Corporation and its affiliates warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at [www.thermofisher.com/us/en/home/global/terms-and-conditions.html](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have questions, contact Life Technologies at [www.thermofisher.com/support](http://www.thermofisher.com/support).



Affymetrix Pte Ltd | 7 Gul Circle #2M-01 | Keppel Logistics Building | Singapore 629563

Revision history: Pub. No. MAN0019853 H00

| Revision | Date             | Description                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                    |
|----------|------------------|----------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| H00      | 7 October 2024   | Removed references to SKUs 905424 and 952433 as they have been discontinued.                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                   |
| G.0      | 16 June 2023     | <ul style="list-style-type: none"><li>• Provided alternative kit option, Axiom™ Consumables Kit for QC (902909), for 903014 and 902801.</li><li>• Updated Part No. 203211, Axygen™ 96-well Round Deepwell Plate, to new Part No. 203367, for the same plate description.</li><li>• Added alternative replacement for the Beckman Coulter™ Biomek™ i7 automated workflow for Axygen™ 384-well Clear 240 µL Polypropylene Deep Well Microplate (non-barcoded), 203103 with Abgene™ 384-Well 250 µL Polypropylene Storage Plate (non-barcoded), 203251.</li><li>• Added information on whether the 384 Deep Well plates are barcoded or non-barcoded in the Beckman Coulter™ automated workflows. Added detailed description of which of these plates are barcoded and non-barcoded in footnote 4 under the Axiom automated workflow for Beckmn Biomek.</li></ul> |
| F.0      | 15 December 2022 | <ul style="list-style-type: none"><li>• Updated several "N/A" fields with the appropriate part number.</li><li>• Deleted PerkinElmer™ Deepwell plate from the PharmacoScan™ NIMBUS® section.</li></ul>                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                         |

The information in this guide is subject to change without notice.

**DISCLAIMER:** TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

**Important Licensing Information:** These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2021-2024 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. Axygen is a trademark of Coming Incorporated. Beckman Coulter and Biomek are trademarks of Beckman Coulter, Inc. Bio-Rad and Hard-Shell are registered trademarks of Bio-Rad Laboratories, Inc. Eppendorf and DeepWell are trademarks of Eppendorf AG. NIMBUS is a trademark of Hamilton Comp. PerkinElmer is a registered trademark of PerkinElmer, Inc. Windows is either a registered trademark or trademark of Microsoft Corporation in the United States and/or other countries.