

Axiom™ NIMBUS™ 2.0 Consumables Kit v2

Catalog Number 952437

Pub. No. MAN0019869 Rev. C.0

Product description

The Axiom™ NIMBUS™ 2.0 Consumables Kit v2 (Cat. No. [952437](#)) contains plastic consumables that are required to run target preparation on the Applied Biosystems™ NIMBUS™ using the Axiom™ 2.0 Reagent Kit (Cat. No. [901758](#)). The kit contains consumables sufficient for 4 × 96 reactions runs. See the appropriate user guide for instructions about using the consumables.

Contents and storage

Component	Part. No. ^[1]	Number		Storage
		Per run	Per kit	
96-well full-skirt plate	203023	7	30	Room temperature
96 half-skirt plate	203244	1	4	
Square deepwell plate	203016	3	12	
Round deepwell plate ^[2]	203367	2	8	
Square 1.2-mL plate	203031	1	4	
4-column reservoir	203056	9	36	

^[1] Plate part numbers are provided for identification purposes only.

^[2] The 96-well round deepwell plate in this kit must be used with the correct thermal shake adapter as it differs from the v1 kit. For details, see the *Axiom™ 2.0 Assay 96-Array Format Automated Workflow User Guide—Applied Biosystems™ NIMBUS™ Target Preparation Instrument* (MAN0017740).

Note: This consumables kit does not include a 96-well UV plate. This plate is in the Axiom™ Consumables Kit for QC (Cat. No. [902909](#)).

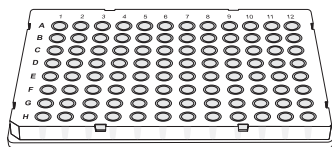


Figure 1 96-well full-skirt plate.

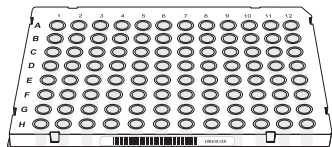


Figure 2 96 half-skirt plate.

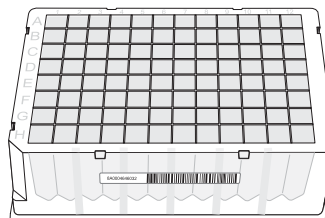


Figure 3 Square deepwell plate.

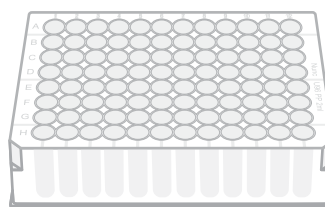


Figure 4 Round deepwell plate.

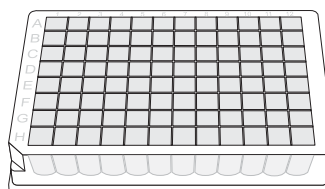


Figure 5 Square 1.2-mL plate.

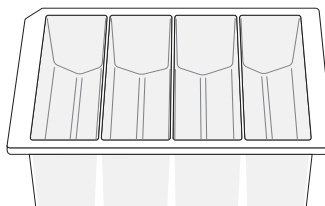


Figure 6 4-column reservoir.

Customer and technical support

Visit [thermofisher.com/support](https://www.thermofisher.com/support) for the latest service and support information.

- Worldwide contact telephone numbers
- Product support information
 - Product FAQs
 - Software, patches, and updates
 - Training for many applications and instruments

- Order and web support
 - Product documentation
 - User guides, manuals, and protocols
 - Certificates of Analysis
 - Safety Data Sheets (SDSs; also known as MSDSs)
- Note:** For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) 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. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



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

Revision history: Pub. No. MAN0019869 C.0

Revision	Date	Description
C.0	25 September 2023	Updated product name.
B.0	10 March 2023	<ul style="list-style-type: none">• Updated 203009 to 203244 and description to "96 half skirt plate".• Updated 203111 to 203367, description to "Round deepwell plate", and quantity to 8.• Updated the respective figure descriptions and images. Updated product name for 902909.

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.

©2023 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.