

Package Insert

CytoScan™ Consumables Kit for NIMBUS™

Intended Use

The CytoScan™ Consumables Kit for NIMBUS™ contains plastic consumables needed to run target preparation on the NIMBUS using the CytoScan™ Assay Kit (P/N 901808). The kit contains consumables sufficient for 4x25 or 2x49 rxn runs. Please refer to the *CytoScan™ Automated Target Preparation Solution on NIMBUS* (P/N 703141) for instructions on using the consumables.

Component	Usage per Run		Number per Kit
	25 Sample	49 Sample	
96-well Semi-skirted Plate	5	7	20
Tube, Screw Cap, No Ribs	6	6	24
1.2 mL Square Well Storage Plate	1	1	4
Reservoir 4 ROW PART PYR PP 73 mL	6	6	25
96-well Full Skirt PCR Plate, Clear	7	7	28
2.2 mL Storage Plate	3	3	12

Storage

The CytoScan™ Consumables Kit for NIMBUS should be stored at room temperature.

Ordering Information

P/N	Product Name	Description
902168	CytoScan™ Consumables Kit for NIMBUS™	Labware sufficient for 4x25 or 2x49 rxn runs

Figure 1. 96-well Semi-skirted Plate

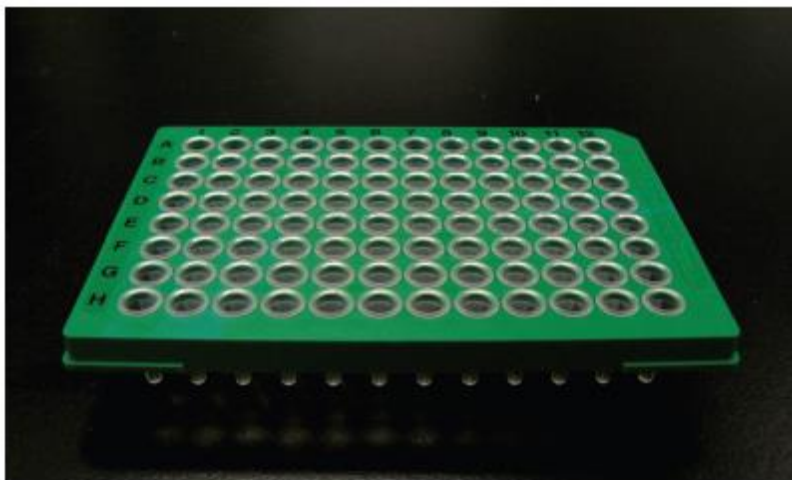


Figure 2. Tube, Screw Cap, No Ribs



Figure 3. 1.2 mL Square Well Storage Plate

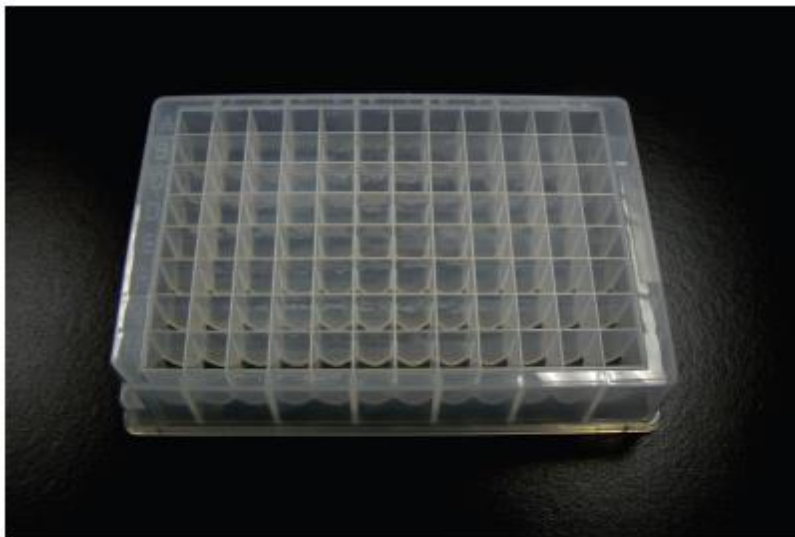


Figure 4. Reservoir 4 ROW PART PYR PP 73 mL



Figure 5. 96-well Full Skirt PCR Plate, Clear

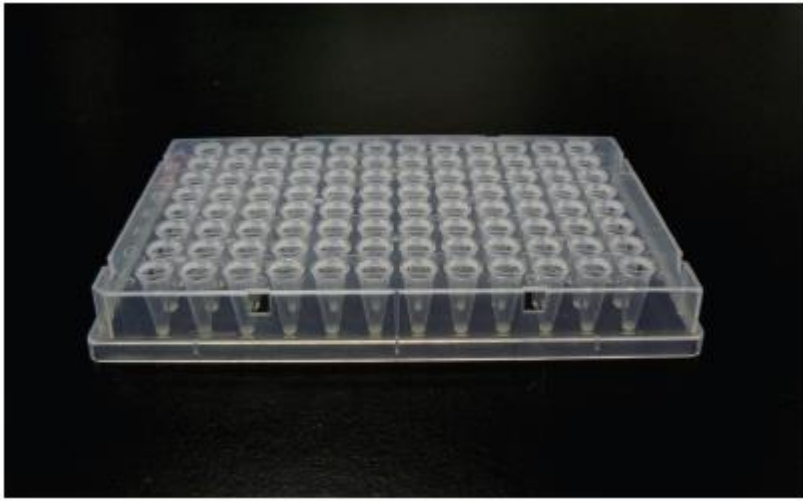


Figure 6. 2.2 mL Storage Plate



Documentation and support

Customer and technical support

Visit thermofisher.com/support for the latest in services and support, including:

- Worldwide contact telephone numbers
- Product support, including:
 - Product FAQs
 - Software, patches, and updates
- Order and web support
- Product documentation, including:
 - 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 found on Life Technologies' web site at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at thermofisher.com/support.

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

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES 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.

Corporate entity: Life Technologies | Carlsbad, CA 92008 USA | Toll Free in USA 1.800.955.6288

©2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. All other trademarks are properties of their respective owners.

P/N 703166

For support visit thermofisher.com/support or email techsupport@lifetech.com

thermofisher.com

23 January 2017

ThermoFisher
SCIENTIFIC