

NEXTFLEX[®] Universal Blockers

(Compatible with Illumina[®] Platforms)

KIT CONTAINS : 16, 48, or 192 RXNS

USER MANUAL FOR :

#NOVA-5143231

#NOVA-5143232

#NOVA-5143233

NEXTFLEX® Universal Blockers

GENERAL INFORMATION _____ 3

 Product Overview 3

 Kit Overview 3

 Warnings and Precautions 3

Usage _____ 4

This product is for research use only.

Not for use in diagnostic procedures.

This manual is proprietary to Revvity, Inc., and intended only for customer use in connection with the product(s) described herein and for no other purpose. This document and its contents shall not be used or distributed for any other purpose without the prior written consent of Revvity. Follow the protocol included with the kit.

Revvity, NEXTFLEX®, NextPrep™, NextPrep-Mag™ are trademarks or registered trademarks of Revvity. All other brands and names contained herein are the property of their respective owners.

GENERAL INFORMATION

Product Overview

The NEXTFLEX® Universal Blockers are designed to prevent concatemerization of library molecules during target capture hybridizations via the adapter sequences that are added to all library molecules before enrichment. Each strand of these adapter sequences can hybridize with the complementary adapter sequence strand of another library molecule during this enrichment hybridization, creating a “daisy-chain” effect between off-target library molecules and the on-target library molecules that are intended targets of the capture probe set used in the enrichment hybridization. By blocking this complementary adapter sequence interaction, the number of on-target reads improves dramatically and the depth of enrichment increases.

The NEXTFLEX® Universal Blockers are designed for use with Illumina® indexed adapters. They are effective in blocking single-indexed or dual-indexed libraries, without being affected by index length or the presence of a unique molecular index (UMI). The NEXTFLEX® Universal Blockers are designed to target both strands of the adapter sequence regions.

The NEXTFLEX® Universal Blockers are compatible with ligation-based workflows, such as the libraries prepared with NEXTFLEX® Unique Dual-Indexed Barcodes. They are also compatible with tagmentation-based workflows, such as Illumina® Nextera™ library prep kits.

Kit Contents, Storage and Shelf Life

The kits contain enough materials to prepare 8, 48, or 96 hybridization reactions. The shelf life is at least 12 months when stored properly. Store at -20°C.

Kit Contents	Cap Color	Amount (8 rxn / 48 rxn / 96 rxn)	Storage Temp.
NEXTFLEX® Universal Blockers	GREY CAP	4.8 / 29 / 58 µL	-20°C

Warnings and Precautions

We strongly recommend that you read the following warnings and precautions. Periodically, optimizations and revisions are made to the components and manual. Therefore, it is important to follow the protocol included with the kit. If you need further assistance, you may contact your local distributor, or contact us at NGS@revvity.com

- Do not use the kit past the expiration date.
- Ensure pipettes are properly calibrated as library preparations are highly sensitive to pipetting error.
- Try to maintain a laboratory temperature of 20°–25°C (68°–77°F).

Version	Date	Description
V21.05	May 2021	Product Launch
V23.11	November 2023	Rebrand to Revvity

Usage

The NEXTFLEX® Universal Blockers can be used with IDT™ xGEN™, Twist Bioscience, and Agilent® SureSelect hybridization capture libraries and workflows.

Recommendations:

IDT™ xGEN™ Hybridization Capture Libraries

Follow Standard protocol of xGEN™ hybridization capture libraries. During hybridization step use 0.6 µL of NEXTFLEX® Universal Blocker per reaction, instead of 2 µL of xGEN™ blocker, independent of library type (Illumina® Truseq™ or Nextera™ type) and index length and/or UMI.

Blocker Master Mix Component	Volume per reaction
Human Cot DNA	5 µL
NEXTFLEX® Universal Blockers	0.6 µL
TOTAL	5.6 µL

Twist Bioscience Hybridization Capture Libraries

Follow Standard protocol of Twist Bioscience hybridization capture libraries. During hybridization step use 0.6 µL of NEXTFLEX® Universal Blocker per reaction, instead of 8 µL of Twist Universal Blocker, independent of library type (Illumina® Truseq™ or Nextera™ type) and index length and/or UMI.

Blocker Master Mix Component	Volume per reaction
Dried Indexed Library Pool	-
Blocker Solution	5 µL
NEXTFLEX® Universal Blockers	0.6 µL
Nuclease-free Water	7.4 µL
TOTAL	13 µL

Agilent® SureSelect Hybridization Capture Libraries

Follow Standard protocol of Agilent® SureSelect hybridization capture libraries. During hybridization step use 0.6 µL of NEXTFLEX® Universal Blocker per reaction, instead of 0.6 µL of Block 3, independent of library type (Illumina® Truseq™ or Nextera™ type) and index length and/or UMI.

Blocker Master Mix Component	Volume per reaction
SureSelect Indexing Block 1	2.5 µL
SureSelect Block 2	2.5 µL
NEXTFLEX® Universal Blockers	0.6 µL
TOTAL	5.6 µL



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.

www.revvity.com

revvity