

Ovation[®] Ultralow Library System V2

www.nugen.com

A single workflow for all of your DNA samples

Ovation[®] Ultralow Library System V2 provides a simple and fast solution for producing libraries that can be used in a broad range of next generation sequencing applications. Using our proprietary **DimerFree** technology allows for efficient library preparation with virtually no adaptor dimers at input levels ranging from 10 pg to 100 ng.

The Ovation Ultralow V2 System is well published across a wide variety of sample types including purified DNA generated from cell lines, fresh and FFPE tissue, liquid biopsy and cell-free DNA. The kit is compatible with most commercially available hybrid capture methods.

Why use Ovation Ultralow Library System V2?

Ultralow V2 offers several unique features:

- 1. Short, simple, automatable workflow:** library construction in as little as 3 hours (2 hours for PCR-free).
- 2. Broad range of input for any sample type:** **DimerFree** technology enables one method to sequence all DNA samples from low to high input, intact or fragmented DNA.
- 3. Accurate variant detection:** 96 Unique Dual Indexes using 192 unique sequences and edit-3 design.
- 4. Simpler workflow:** no hidden sample normalization or adaptor dilution steps required.
- 5. No adaptor dimers:** get more informative reads compared to competitors without multiple bead clean-ups.
- 6. Obtain broad coverage:** high fidelity amplification across a broad range of GC content.

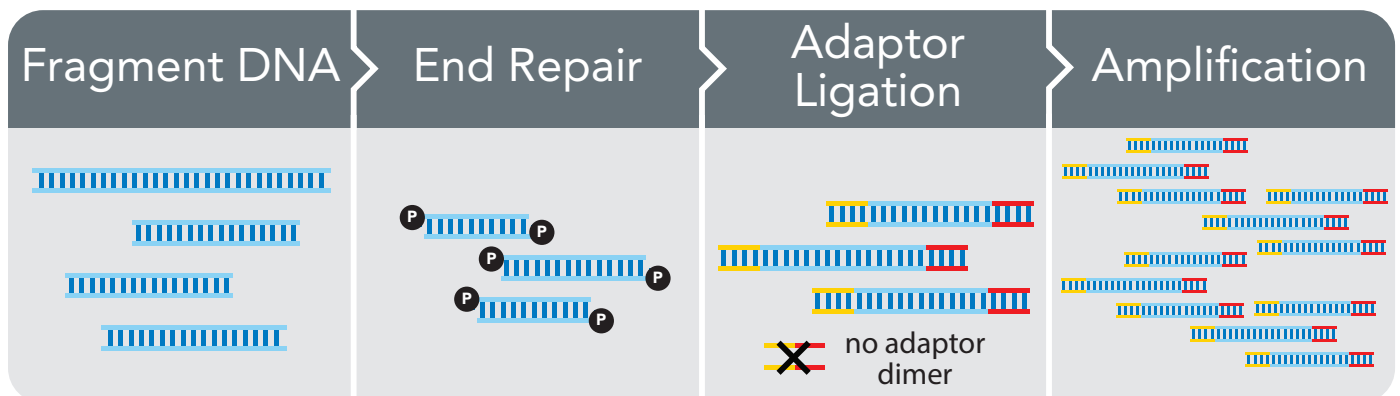


Figure 1. Ovation Ultralow System V2 is a simple, fast and scalable workflow that produces clean libraries with no adaptor dimers.

Applications and Sample Types

- Whole genome sequencing
- Exome sequencing
- ChIP-Seq
- Microbiome
- Liquid biopsy
- FFPE DNA
- cDNA

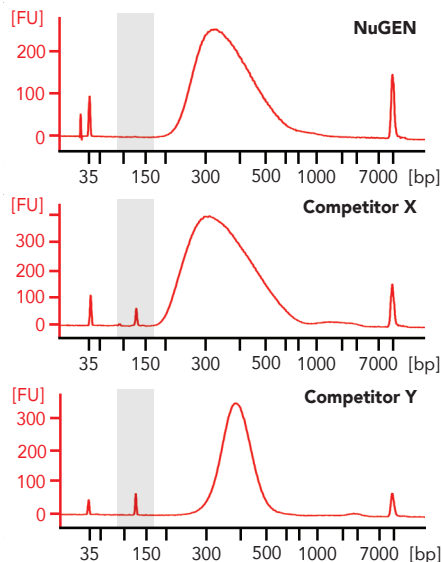


Figure 2. No detectable adaptor dimers. 1 ng gDNA created from Ultralow System V2 vs. Competitors X and Y analyzed on Agilent Bioanalyzer. Adaptor dimers are highlighted in the grey box.

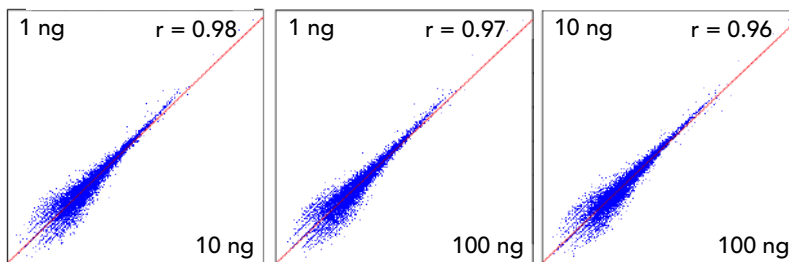


Figure 3. High concordance across a broad input range without the need to adjust adaptor concentrations. Scatterplots of the log FPKM and corresponding pearson R values are shown for Ultralow Library Systems V2 libraries constructed from 1, 10 or 100 ng of double-stranded cDNA made from Human Brain reference RNA.

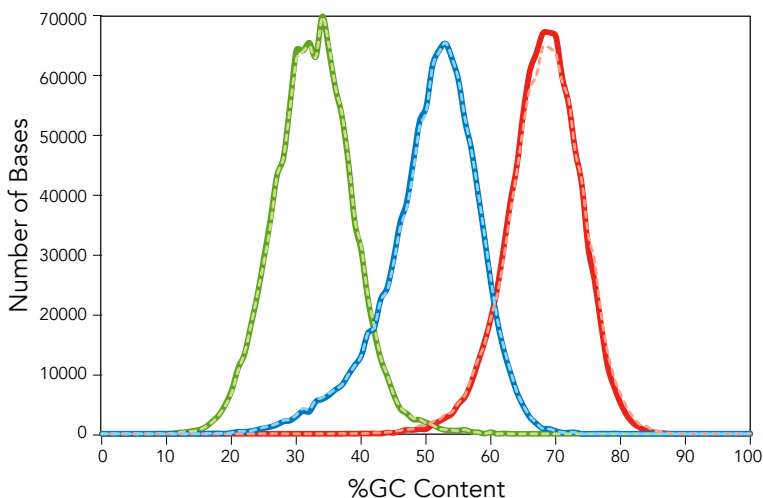


Figure 4. Even read distribution across a broad GC range. 1 ng of gDNA from *E. coli* (blue, 51% GC), *R. sphaeroides* (red, 69% GC) or *S. aureus* (green, 33% GC) were sequenced with Ultralow Library Systems V2. Experimental GC distribution (solid lines) closely matches theoretical distribution (dashed lines) from a GC of less than 20% to greater than 80%.

Ordering Information	No. of Reactions	Part No.
Ovation Ultralow System V2	8	0344NB-08
Ovation Ultralow System V2, includes purification beads	32	0344-32
Ovation Ultralow System V2	32	0344NB-32
Ovation Ultralow System V2 1-96	96	0344NB-A01
Ovation Ultralow System V2 with Unique Dual Indexes	96	9149-A01



www.nugen.com



NuGEN Technologies, Inc.

Headquarters USA

201 Industrial Road, Suite 310
San Carlos, CA 94070 USA
Toll Free Tel: 888.654.6544
Toll Free Fax: 888.296.6544
custserv@nugen.com
techserv@nugen.com

Europe

P.O. Box 109,
9350 AC Leek
The Netherlands
Tel: +31-13-5780215
Fax: +31-13-5780216
europe@nugen.com

Worldwide

For our international distributors contact information, visit our website.
worldwide@nugen.com



©2017-2018 NuGEN Technologies, Inc. All rights reserved. The Encore®, Ovation® and Applause® families of Products and methods of their use are covered by several issued U.S. and International patents and pending applications (www.nugen.com). NuGEN, Allegro, Celero, NuQuant, SoLo, Metaplex, Ovation, SPIA, Ribo-SPIA, Applause, Encore and Imagine More From Less are trademarks or registered trademarks of NuGEN Technologies, Inc.

Other marks appearing in these materials are marks of their respective owners.

For research use only.

M01468 v1.1