

PRODUCT INFORMATION Thermo Scientific Verso cDNA Synthesis Kit

#AB-1453/A

40 rxns

Lot _

Expiry Date _

Ordering Information

	#AB-1453/A 40 rxns of 20 µL	#AB-1453/B 100 rxns of 20 μL
Verso Enzyme Mix	40 μL	100 μL
5X cDNA Synthesis Buffer	500 μL	500 μL
Anchored Oligo dT (500 ng/µL)	40 μL	100 μL
Random Hexamer (400 ng/µL)	40 μL	100 μL
dNTP Mix (5 mM each)	200 μL	200 μL
RT Enhancer	40 μL	100 μL

Store at -20 °C



Description

Thermo Scientific Verso cDNA Kit supplies all the reagents to generate high yields of full-length cDNA from all RNA types.

Verso™ Enzyme Mix includes Verso Reverse
Transcriptase, which is active at high temperatures, is
highly sensitive and can generate long cDNA strands. This
mix also contains RNase inhibitor to protect RNA
templates from degradation.

5X cDNA Synthesis Buffer, a proprietary reaction buffer which has been optimized to improve reverse transcription across a wide range of templates.

Anchored Oligo dT primers and **Random Hexamers** provide flexible RNA priming methods for cDNA synthesis.

RT Enhancer is included to remove contaminating DNA, eliminating the need for DNAse I treatment. It degrades double stranded DNA during the transcription of RNA and is inactivated after 2 minutes at 95 °C.

Verso Reverse Transcriptase

Verso is an RNA-dependent DNA polymerase with a significantly attenuated RNase H activity. Verso can synthesize long cDNA strands, up to 11 kb, at a temperature range of 42 °C to 57 °C. The recommended amount of total RNA to use is between 1 pg and 1 µg.

Storage Conditions

Store at -20 °C until ready for use. Avoid repeated freeze thawing.

Additional Info

The use of disposable gloves, RNase and DNase free filter tips and plastics is recommended.

If DNase I treatment has been performed, RT Enhancer is not required.

Tips before use

Thaw the reagents on ice. Mix and spin down the solutions before use to recover the maximum amount. **Do not vortex the Verso Enzyme Mix.**

Briefly centrifuge to avoid bubbles within the wells. Always include a no template control (NTC) and a no enzyme control (NEC).

Protocol

Example of reaction mix preparation.

The volume of each component is for a 20 µL final reaction.

	Volume	Final Concentration
5X cDNA synthesis buffer	4 µL	1X
dNTP Mix	2 µL	500 µM each
RNA Primer*	1 µL	
RT Enhancer	1 µL	
Verso Enzyme Mix	1 µL	
Template (RNA)	1-5 µL	1 ng
Water, nuclease-free (#R0581)	To 20 μL	
Total volume	20 µL	

^{*}It is recommended that RNA primers be added to the <u>final</u> 1X reaction as follows: 1 μ L of anchored oligo dT (orange cap) <u>or</u> 1 μ L of random hexamers (blue cap) <u>or</u> 1 μ L of a blend of random hexamers and anchored oligo-dT 3:1 (v/v) <u>or</u> gene-specific primer (to final concentration of 0.5 – 2 μ M).

Anchored oligo dT is not suitable for use with most prokaryotic RNA. In these cases, random hexamers or gene-specific primers are recommended.

Example of reverse transcription cycling program:

	Temp.	Time	Number of cycles
cDNA synthesis*	42 °C	30 min	1 cycle
Inactivation	95 °C	2 min	1 cycle

^{*}Depending on the length of template and degree of secondary structure, the efficiency of the first strand synthesis may be improved by optimizing temperature and time (42-57 °C for 5-60 minutes).

CERTIFICATE OF ANALYSIS

Verso cDNA Kit is tested functionally for use in RT-PCR.

Quality authorized by:

Jurgita Zilinskiene

LIMITED USE LABEL LICENSE: Internal Research and Development Use Only.

The purchase of this product conveys to the buyer the limited, non-exclusive, non-transferable right (without the right to resell, repackage, or further sublicense) to use this product for internal research and development purposes. No other license is granted to the buyer whether expressly, by implication, by estoppel or otherwise. In particular, the purchase of the product does not include or carry any right or license to use, develop, or otherwise exploit this product commercially and no rights are conveyed to the buyer to use the product or components of the product for purposes including but not limited to provision of services to a third party, generation of commercial databases or clinical diagnostics. This product is sold pursuant to authorization from Thermo Fisher Scientific and Thermo Fisher Scientific reserves all other rights. For information on purchasing a license for uses other than internal research and development purposes, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies Inc., 5791 Van Allen Way, Carlsbad, California 92008.

PRODUCT USE LIMITATION

This product is developed, designed and sold exclusively for research purposes and in vitro use only. The product was not tested for use in diagnostics or for drug development, nor is it suitable for administration to humans or animals. Please refer to www.thermoscientific.com/onebio for Material Safety Data Sheet of the product.

© 2015 Thermo Fisher Scientific, Inc. All rights reserved. All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.