



PRODUCT INFORMATION

Random Hexamer Primer

5' -d (NNNNNN) -3' N = G, A, T or C

Pub. No. MAN0013111

Rev. Date 05 October 2017 (Rev. B.00)

#SO142

Assembling Lot 00000000

Filling Lot 00000000

Expiry Date MM.YYYY

Store at -20 °C

Components	#SO142
Random Hexamer Primer, 100 µM (0.2 µg/µL)	120 µL (24 µg)

www.thermofisher.com

For Research Use Only. Not for use in diagnostic procedures.

Description

The Random Hexamer Primer is a mixture of single-stranded random hexanucleotides with 5'- and 3'-hydroxyl ends.

The primer is supplied as a ready-to-use, 20X concentrated aqueous solution.

Application

First strand cDNA synthesis.

CERTIFICATE OF ANALYSIS

Functionally tested in RT-PCR using 100 fg of control GAPDH RNA.

Quality authorized by:

Jurgita Zilinskiene

Rev.6



53

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., 5781 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.thermofisher.com for Material Safety Data Sheet of the product.

© 2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.