

## PRODUCT INFORMATION

# Anchored Oligo dT Primers

#AB-1247      100 µL

Lot \_      Exp. Date \_

Store at -20°C



[www.thermoscientific.com/onebio](http://www.thermoscientific.com/onebio)

## Description

Anchored Oligo-dT Primers are specifically designed for synthesising long cDNAs from mRNAs carrying poly-A tails. Since this primer only anneals to the poly-A tail junction of the mRNA template, it eliminates transcription through the poly-A tail providing more effective cDNA synthesis.

**Concentration:** 500 ng/µL

## 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

## 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](http://www.thermoscientific.com/onebio) for Material Safety Data Sheet of the product.

© 2012 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

