

TaqMan® Copy Number Reference Assays

Insert PN 4403453 Rev. D

For research use only. Not intended for human or animal therapeutic or diagnostic use.

These reference assays are optimized for use with the TaqMan® Copy Number Assays to detect and measure copy number variation in the human or mouse genome. Refer to the *TaqMan® Copy Number Assays Protocol* (PN 4397425) and the *CopyCaller™ Software User Guide* (PN 4400042) for detailed information on the experimental procedure. The reference assays are species-specific.

Table 1 Human assays

| Part number | Part | Volume | Number of reactions (20 µL reaction volume) | Concentration | Storage conditions |
|-------------|---|--------------------------|---|---------------|---|
| 4403326 | TaqMan® Copy Number Reference Assay RNase P <ul style="list-style-type: none"> Two unlabeled PCR primers TAMRA™ probe (VIC® dye-labeled) | 1 tube (750 µL) | 750 reactions | 20X | Store at – 15 °C to – 25 °C. Protect from light. Minimize freeze-thaw cycles. |
| 4403328 | | 4 tubes (750 µL each) | 3000 reactions | | |
| 4403316 | TaqMan® Copy Number Reference Assay TERT <ul style="list-style-type: none"> Two unlabeled PCR primers TAMRA™ probe (VIC® dye-labeled) | 1 tube (750 µL) | 750 reactions | | |
| 4403315 | | 4 tubes (750 µL each) | 3000 reactions | | |

Table 2 Mouse assays

| Part number | Part | Volume | Number of reactions (20 µL reaction volume) | Concentration | Storage conditions |
|-------------|--|--------------------------|---|---------------|---|
| 4458366 | TaqMan® Copy Number Reference Assay, Mouse, Tfrc <ul style="list-style-type: none"> Two unlabeled PCR primers TAMRA™ probe (VIC® dye-labeled) | 1 tube (750 µL) | 750 reactions | 20X | Store at – 15 °C to – 25 °C. Protect from light. Minimize freeze-thaw cycles. |
| 4458367 | | 4 tubes (750 µL each) | 3000 reactions | | |
| 4458368 | TaqMan® Copy Number Reference Assay, Mouse, Tert <ul style="list-style-type: none"> Two unlabeled PCR primers TAMRA™ probe (VIC® dye-labeled) | 1 tube (750 µL) | 750 reactions | | |
| 4458369 | | 4 tubes (750 µL each) | 3000 reactions | | |

Note: For safety and biohazard guidelines, refer to the “Safety” section in the *TaqMan® Copy Number Assays Protocol* (PN 4397425). Read the Safety Data Sheet (SDS) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

For research use only. Not intended for human or animal therapeutic or diagnostic use.

NOTICE TO PURCHASER: PLEASE REFER TO THE TAQMAN COPY NUMBER ASSAYS® PRODUCT INSERT AND PROTOCOL FOR LIMITED LABEL LICENSE OR DISCLAIMER INFORMATION.

The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners. AmpliTaq Gold and TaqMan are registered trademarks of Roche Molecular Systems, Inc.

© Copyright 2010, Life Technologies Corporation. All rights reserved.

Part Number 4403453 Rev. D 10/2010



Headquarters

5791 Van Allen Way | Carlsbad, CA 92008 USA
Phone 760.603.7200
www.lifetechnologies.com

International Sales

For our office locations please call the division headquarters or refer to our Web site at
www.appliedbiosystems.com/about/offices.cfm