

LANCE® *Ultra***Human c-Myc Detection Kit****Product number:** TRF1111C **Lot Number:** 3310354**Material provided:****Kit Format:** TRF1111C: 500 assay points TRF1111M: 10 000 assay points

The number of assay points is based on an assay volume of 20 µL in 384-well assay plates using kit components at the recommended assay concentrations.

Standard Format: TRF1111S: 100 µL at 10mM**Manufacturing date:** 7/02/2024 **Document version:** 1**Product Information**

- Application:** This kit is designed for the quantitative determination of c-Myc in media using a homogeneous LANCE *Ultra* assay (no wash steps).
- Kit contents:** The kit contains 5 components: LANCE *Ultra* Eu-labeled Streptavidin, LANCE *Ultra* ULight-labeled Anti-c-Myc Antibody stored in TSA, 0.1% BSA, c-Myc Analyte, Biotinylated c-Myc Tracer, and 5X *Ultra* HiBlock Buffer.
- Sensitivity:** Lower Detection Limit (LDL): 46.1 nM
Lower Limit of Quantification (LLOQ): 235.8 nM
EC₅₀: 1 346.4 nM
- Dynamic Range:** 46.1– 300,000 nM
- Storage:** Store kit in the dark at +4°C. Store analyte at -20°C upon receipt. Store reconstituted analyte at -20°C.
- Stability:** This kit is stable for at least 12 months from the manufacturing date when stored in its original packaging and the recommended storage conditions.

Quality Control

Lot to lot consistency is confirmed in an LANCE *Ultra* assay. EC50 and LDL were measured on the, EnVision Multilabel Plate Reader equipped with TR-FRET option and a TRF laser using the protocol described in this technical data sheet. We certify that these results meet our quality release criteria. Assay performance may vary between lots and by the instrument used.

EC50: 1.648 µM
LDL: 106.142 nM
LLOQ: 619.413 nM
Min Counts: 958
Max Counts: 12,387

The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.

TRF1111-R Rev01

The Revvity logo is displayed in a lowercase, sans-serif font.

Revvity, Inc.
940 Winter Street
Waltham, MA 02451 USA

(800) 762-4000 www.revvity.com

For a complete listing of our global offices, visit www.revvity.com
Copyright ©2023, Revvity, Inc. All rights reserved.