

LANCE® *Ultra***EGFR Positive Control Lysate****Product number:** TRF4015S**Lot Number:** 3254296**Kit Format:** TRF4015S: 500µL frozen cell lysate**Product Information****Application:** LANCE *Ultra* EGFR Positive Control Lysate. This product is designed to be used in combination with LANCE *Ultra* Cellular Detection kits: TRF4015**Formulation:** A431 cells treated with EGF for 10 minutes prior to lysis with LANCE *Ultra* Lysis Buffer 1 at room temperature for 30 minutes and then frozen.**Storage:** Store LANCE *Ultra* Lysate at -80 °C**Stability:** This product is stable for at least 3 months from the date of receipt when stored in its original packaging and the recommended storage conditions.

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.

TRF4015S-R Rev01