

LANCE® *Ultra*

## eIF2A (S51) Cellular Detection Kit

**Product number:** TRF4008C      **Lot Number:** 3220348

**Kit Format:** TRF4008C: 500 assay points      TRF4008M: 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.

**Manufacturing date:** November 7, 2023      **Document version:** 1

### Product Information

**Application:** This kit is designed for the detection of phosphorylated eIF2A (s51) in cell lysates using a homogeneous LANCE *Ultra* assay (no wash steps).

**Kit contents:** The kit contains 4 components: LANCE *Ultra* Eu-labeled Anti-eIF2A (s51) Antibody stored in TSA, 0.1% BSA, LANCE *Ultra* ULight-labeled Anti-eIF2A Antibody stored in TSA, 0.1% BSA, 10X LANCE Detection Buffer, and 10X LANCE *Ultra* Lysis Buffer 2.

**Typical Performance:** Signal/Background: 25      Z': 0.96

**Storage:** Store kit in the dark at +4°C.

**Stability:** This kit is stable for at least 6 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. S/B and Z' 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.

SB: 17.5  
Z': 0.871

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.

TRF4008-R Rev01