

Alpha SureFire® Ultra™ Multiplex

p-Rb (Ser780) + Total Rb

Part numbers: MPSU-PTRB-K-L

Lot Number: A1664

Manufacturing date: December 02, 2024



CONTENTS:

p-Rb (Ser780) + Total Rb
Positive Control (lyophilised)

Batch #:

N1409

STORAGE / HANDLING

Positive Control Lysate store at -20°C. Once reconstituted, lysate can generally be stored effectively at -20°C in single use aliquots but should be tested on a case by case basis.

KIT STABILITY:

Stable for at least 4 months from the manufacturing date when stored under recommended storage conditions.

CONTROL LYSATE EXPECTED SIGNAL:

Reconstituted control lysate diluted to 6.25 % should give a signal window over buffer alone of at least the following:

p-Rb (Ser780)	170 - fold	Total Rb	40 - fold
----------------------	-------------------	-----------------	------------------

This recommended dilution falls within the linear range of the assay. The control lysate signal window can be influenced by lysate solubility, differing lab conditions, reader settings and reader performance.

This product is not for resale or distribution except by authorized distributors.

For more information on protocols, please refer to the generic AlphaLISA SureFire Ultra Assay Kits Manual, available from www.revvity.com

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.