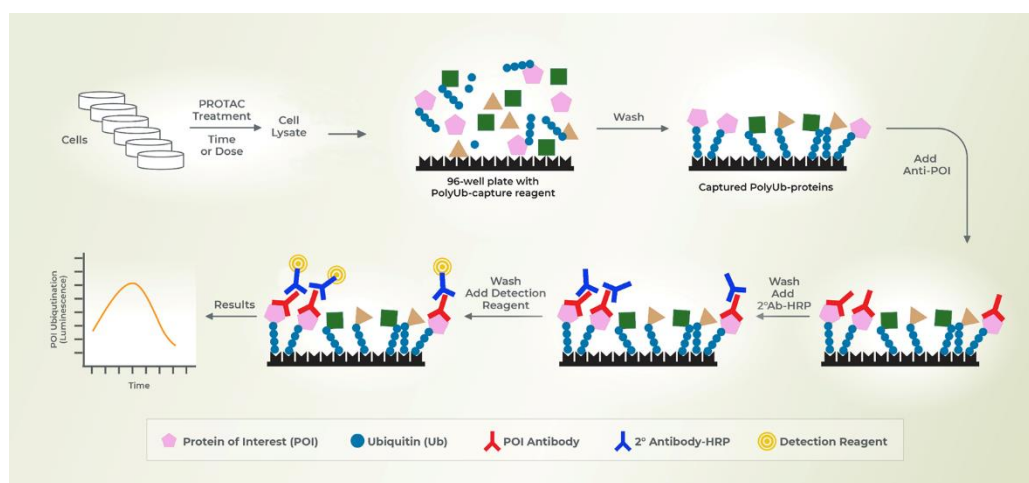


K63 Ubiquitin Linkage ELISA Kit (Chain Selective)

MANUAL

K63 Ubiquitin Linkage ELISA Kit (Chain Selective) Catalog Number: PA630



all products are for research use only • not intended for human or animal diagnostic or therapeutic uses
LifeSensors, Inc., 271 Great Valley Parkway, Malvern PA 19355 • (p) 610.644.8845 (f) 610.644.8616
techsupport@lifesensors.com • www.lifesensors.com • sales@lifesensors.com
Copyright © 2025 LifeSensors, Inc. All Rights Reserved

K63 Ubiquitin Linkage ELISA Kit (Chain Selective)

Cat. # PA630

BACKGROUND

The Ubiquitin Proteasome System (UPS) is a dynamic process involved in various cellular functions, including cell signaling, DNA repair, and proteasomal degradation. Ubiquitin can form diverse chains on its seven lysine residues, with K48 and K63-linked polyubiquitination being a crucial post-translational modification. While K48 polyubiquitination predominantly drives protein degradation and maintains cellular homeostasis, K63 linked polyubiquitination plays pivotal role in various cellular processes, primarily associated with trafficking, signaling pathways, DNA repair mechanisms and the regulation of protein-protein interactions.

INTENDED USES

LifeSensors' K63 Ubiquitin Linkage ELISA kit is intended to enrich and quantitate K63 ubiquitination in cellular and tissue lysates or in in vitro ubiquitination reactions. This kit enables relative and absolute quantitation of K63 ubiquitination on target proteins in cellular and tissue lysates under various treatments (PROTACs, Molecular Glues, MG132, etc) and conditions. The plate is designed for research use only and is not intended for human or animal diagnostic or therapeutic applications.

BENEFITS

- Replaces labor-intensive, semi-quantitative Western blot methods for examining K63 polyubiquitination and protein degradation.
- Offers quantitative, reproducible results.
- Facilitates high-throughput screening for compound library processing and structure-activity relationship (SAR) studies.

PRINCIPLE OF THE ASSAY

Microtiter plate strips, pre-coated with a proprietary K63-TUBE reagent (assay plate) are used to capture K63-linked polyubiquitinated proteins in cell or tissue lysates or during an in vitro ubiquitination reaction. Non-polyubiquitinated proteins or those ubiquitinated on lysines residues other than K63 are either excluded from binding or removed by washing. The K63-ubiquitinated target Protein of Interest (POI) is detected with primary antibody against the POI (not included) and a secondary antibody conjugated with-HRP (included) allowing for detection by chemiluminescence (Figure 1). Thus, the signal generated in this "sandwich" ELISA assay is a quantitative measure of K63-linked ubiquitination of target proteins.

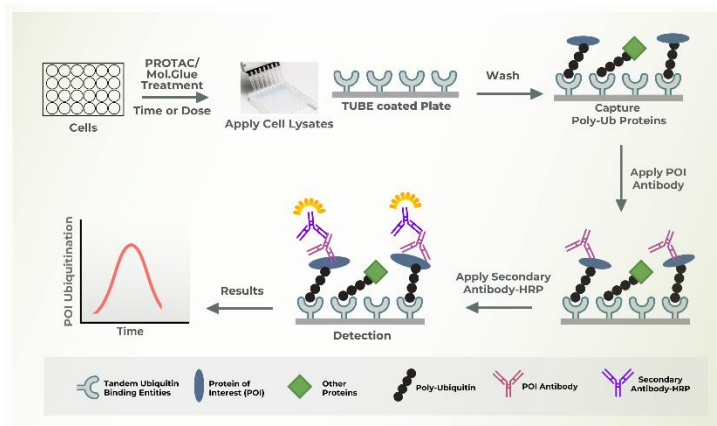


Figure 1. Schematic of K63 Ubiquitin Linkage ELISA Assay.

K63 Ubiquitin Linkage ELISA Kit (Chain Selective)

COMPONENTS INCLUDED IN THE KIT

1. **K63-Ubiquitin Chain Assay 96-well strip well plate**
Note: Do not thaw the assay plate rapidly to room temperature. We recommend placing the assay plate at 4°C for 1 hour prior to transferring to room temperature.
2. **Blocking Concentrate (BC, 5X) –1 x 12 mL**
3. **Detection Reagent1 (DR1, 2X)- 1x 1.0 mL of Hydrogen Peroxide**
4. **Detection Reagent 2 (DR2, 2X)–1 x 1.0 mL of Luminol**
5. **UPS Inhibitor cocktail (10X)- 1x 1.0 mL (For adding to the Lysis Buffer)**
6. **Proteasome Inhibitor (10X)- 1x 1.0 mL (For adding to the Lysis Buffer)**
7. **Plastic plate seals –2 x**

Positive Control Reagents:

8. **Biotin-K63-Poly-Ubiquitin Chain (20X): - 1 x 25 µL**
9. **Streptavidin HRP (100X): 1 x 10 µL**

STORAGE AND STABILITY

Store Coated Plate at -80°C. Store Blocking Concentrate and Detection Reagents at 4°C. Store the UPS and Proteasome Inhibitor Cocktails at -20°C. Plate seal can be stored at room temperature. Avoid multiple freeze/thaw cycle. All components are stable for at least 6 months if stored at the proper temperature.

ADDITIONAL ITEMS REQUIRED BUT NOT INCLUDED IN THE KIT

1. **Purified Target Protein of Interest if performing in vitro ubiquitination**
2. **PROTACs, Molecular Glues, MG132 etc for cell treatment or for in vitro ubiquitination reaction**
3. **Anti-POI Specific Primary Antibodies (ELISA compatible and highly specific)**
4. **WASH BUFFERS: 1X Phosphate Buffered Saline, 0.1% Tween (PBS-T)**
5. **Phosphatase inhibitors if studying phosphorylated target proteins of interest**
6. **LUMINESCENCE CAPABLE PLATE READER**
7. **15 mL centrifuge tubes**
8. **Polypropylene plate for preparing compound dilutions (Optional)**
9. **Multi-channel Pipettors and Automatic Plate Washer (Optional)**
10. **Disposable gloves, pipette tips and reagent reservoirs**

OPTIONAL

Decomplexing Buffer - SKU: UE-1003-2000. This buffer is available for purchase on the LifeSensors website [here](#).

For some targets, the decomplexing agent enhances the signal to background. We recommend the use of the decomplexing agent prior to analysis with PA950, PA480 & PA630. The use of the decomplexing agent disrupts any native protein complexes that might be part of ubiquitin complexes. This urea based decomplexing agent results in reduced background signal, resulting in a better signal-to-background ratio.

K63 Ubiquitin Linkage ELISA Kit (Chain Selective)

DETAILED PROTOCOL

Note: It is essential that the entire manual is thoroughly reviewed before starting the procedure.

1. Remove the plate from freezer and reagents from the refrigerator, and place them at 4°C for at least 1 hour. Then allow them to reach room temperature (RT). Incubations are performed at either 4°C or RT (22°-27°C). To avoid cross contamination, do not re-use plate sealers.
2. Prepare cell lysate (described below in Cell Lysis Protocol) and decomplex the lysate with Decomplexing Buffer for 15 min with vortexing at 5 min interval. Add 15-25 µg/well of decomplexed lysate diluted with PBS (Total Volume = 100 µl/well). Seal the plate.
3. Incubate plate for 1.5 - 2 hours at RT with gentle orbital shaking (~300rpm).
4. Wash the assay plate 4 times with PBS-T using 200µL per well, tap out (upside down) excess PBS-T from wells between washes on paper towels or other blotting paper. **NOTE: DO NOT ALLOW WELLS TO DRY COMPLETELY.**
5. Prepare sufficient volume of 1X Blocking Concentrate by diluting 5X stock into 1X PBST.
6. Optional step: If using the supplied positive control (Biotin-K63-polyubiquitin chain) to verify plate performance, add 17.5 µL of the 20X Biotin-K63-polyubiquitin chain into 332.5 µL 1X BC
7. Dilute the anti-POI primary antibody according to antibody manufacturer instructions in (1X) BC.
 - i. *Note: Primary antibody dilution depends on primary antibody efficiency and manufacturer's recommendation. A good starting dilution is 1:1000 or 0.5-1 µg/mL.*
8. Pipette 100 µL of diluted **anti-POI primary antibody** into each well.
9. Optional step: Pipette 100 µL of diluted Biotin-polyubiquitin chain (positive control) into three wells in assay plate. Pipette 100 µL 1X BC into three wells as blank control.
10. Seal the plate and incubate the assay plate for **60 minutes** at **room temperature** with shaking at 300RPM on an orbital shaker.
11. Remove the plate sealer and decant the reagents from the plate. Wash the assay plate 4 times with PBS-T using 200µL per well, tap out (upside down) excess PBS-T from wells between washes on paper towels or other blotting paper. **NOTE: DO NOT ALLOW WELLS TO DRY COMPLETELY.**
12. Dilute HRP-conjugated secondary antibody in 1x Blocking Concentrate and Dilute the Streptavidin-HRP to 1X in 1X BC for positive control detection.
13. Add the diluted secondary Antibody-HRP conjugate 100 µL/well. Optional step: Pipette 100µL of Streptavidin-HRP in 1X BC into wells containing positive control and blank. Incubate plate for 45 minutes at RT with gentle orbital shaking (~300rpm).
 - i. *Note: Secondary Antibody-HRP conjugate should be diluted as recommended by manufacturer. A good starting dilution range is 1:1000 to 1:5000.*
14. Remove the plate sealer and decant the reagents from the plate. Wash the assay plate 4 times with PBS-T using 200µL per well, tap out excess PBS-T from wells between washes.
15. Just before use, mix 800 µL of DR1 and 800 µL of DR2 into 10 mL of ultrapure water (deionized or distilled). Add 100 µL of this detection reagent solution to each well and read immediately with a plate reader optimized for chemiluminescence for 5-10 reads with 1 min intervals.
 - a. **Important Note:** DR1 and DR2 amounts needed will vary based on overall signal and plate reader being used. Optimize the amount of DR1 and DR2 needed by reading with a 1:5 dilution of each reagent and then titrating up (1:10 to 1:20) based on initial raw values.

K63 Ubiquitin Linkage ELISA Kit (Chain Selective)

CELL LYSIS PROTOCOL

1. Aspirate medium completely and rinse cells with ice cold 1X PBS. For adherent cells, add lysis buffer directly to the plate or wells and scrape cells into the lysis buffer (~75µl/well for a six-well plate). For suspension cells, spin down cells at ~1500 rpm for 3-5 minutes at 4°C. Gently aspirate media, wash the pellet with ice cold 1X PBS. Freeze cell pellet at -80°C for long term storage or proceed with lysis on ice.
2. For cell pellet, add RIPA lysis buffer (5-10 times the volume of the pellet, for example, for 100 µL of pellet add 500-1000 µL lysis buffer).
3. Vortex the cell lysate intermittently for 10-15 minutes while keeping the sample on ice to allow efficient lysis.
4. Centrifuge at 13,000 x g for 15-30 min at 4°C.
5. Collect the supernatant (whole cell lysate) and determine protein concentration using either Bradford or BCA method following manufacturer's recommendations.

RIPA Lysis Buffer

50 mM Tris-HCl, pH 7.5

150 mM NaCl

1% NP40

1% Sodium deoxycholate

2 mM EDTA

1 mM PMSF

10% Glycerol

1X UPS inhibitor cocktail (Add just before use)

1X Proteome inhibitor cocktail (Add just before use)

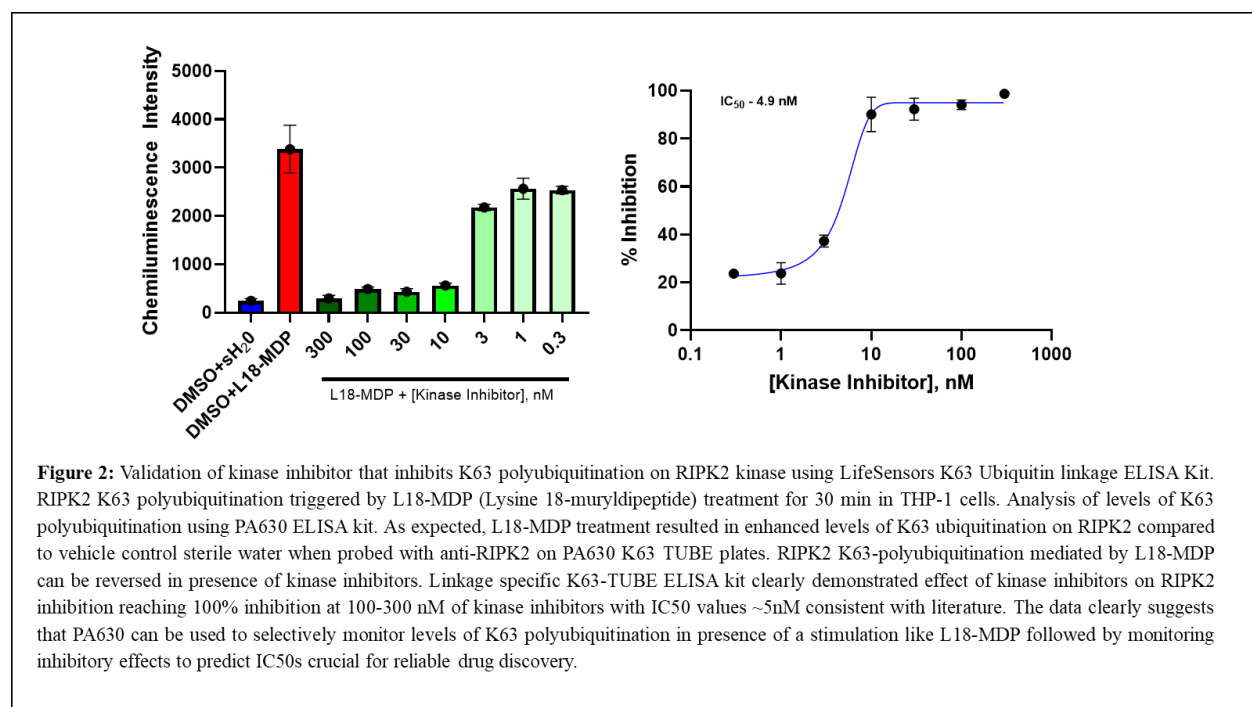
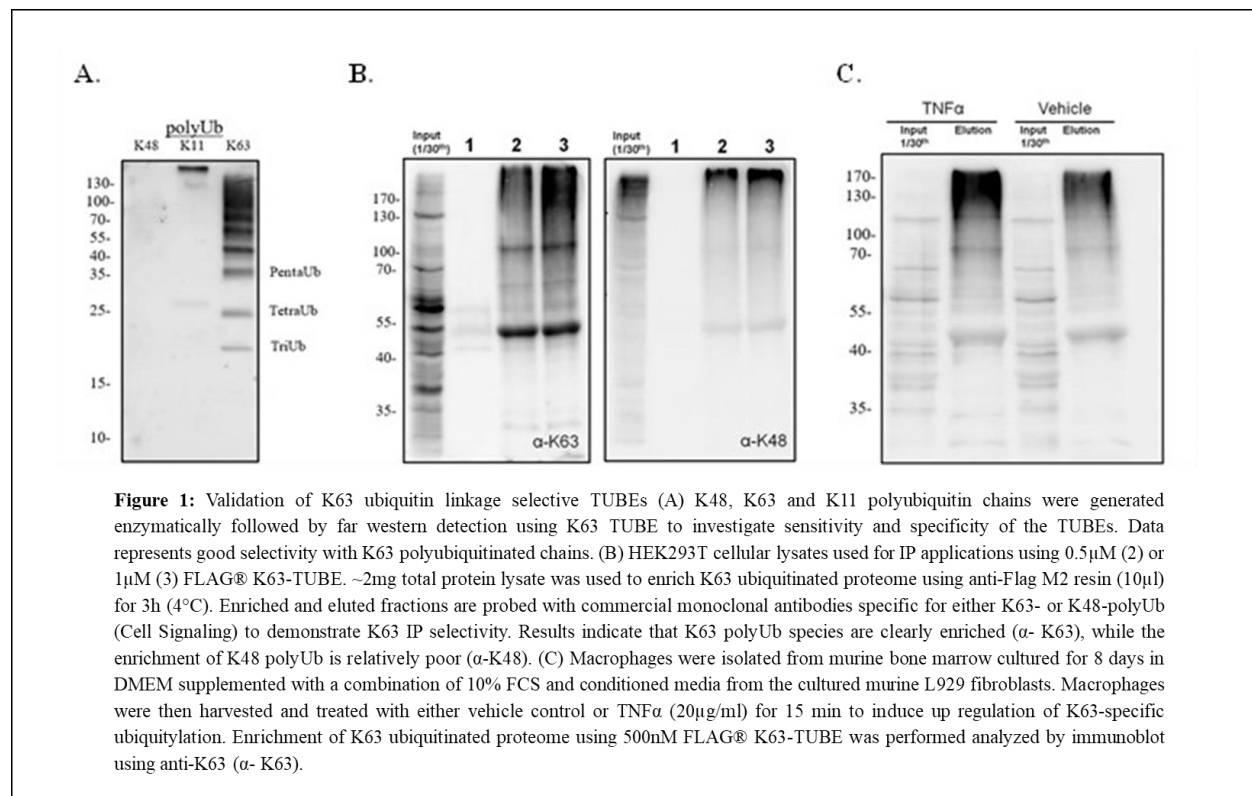
1:500 Protease Inhibitor cocktail (Sigma Cat# P8849) (Add just before use)

NOTES

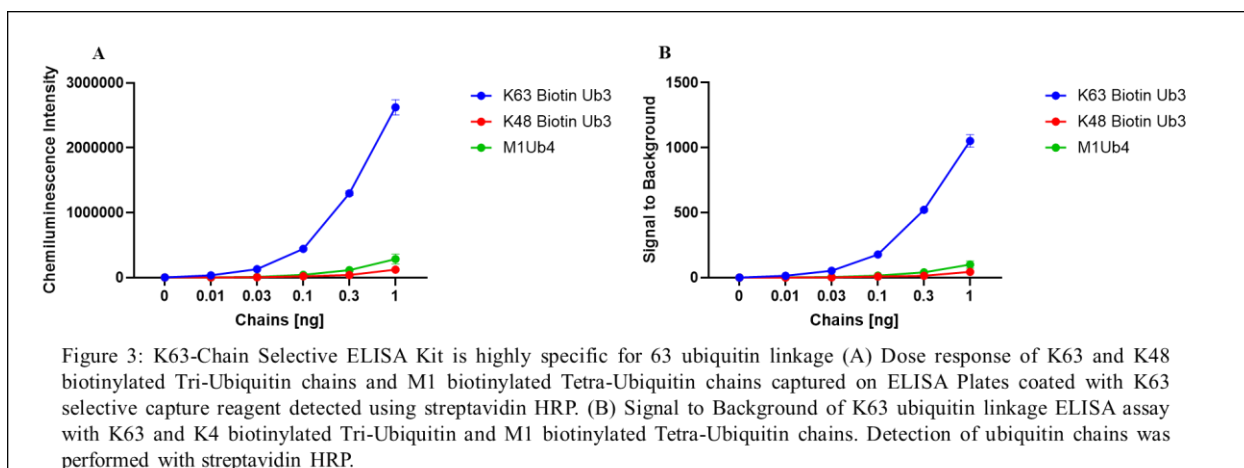
1. For Targeted protein degradation applications - to monitor K63 polyubiquitination of target protein followed by its degradation, please select times before and after D_{max} of PROTAC.
 2. Optimize antibody dilution and specificity of target protein by western blotting. Selection of a good, clean target-specific antibody suitable for sandwich ELISA is essential for the success of this assay.
 3. Optimize lysate concentration depending on target abundance.
 4. See suggested lysis buffer below to have the best polyubiquitination profile of target protein.
 5. Include appropriate controls for best results and interpretation.
-

K63 Ubiquitin Linkage ELISA Kit (Chain Selective)

REPRESENTATIVE DATA



K63 Ubiquitin Linkage ELISA Kit (Chain Selective)



K63 Ubiquitin Linkage ELISA Kit (Chain Selective)

TROUBLESHOOTING GUIDE

Below are common problems, possible causes, and recommended solutions to help you achieve reliable and reproducible results.

1. Weak or No Chemiluminescence Signal

Possible Causes & Solutions:

- Poor primary antibody performance → This can be a major hurdle in getting optimal results. Verify that the 1° antibody is validated for ELISA and titrate as needed. A poor antibody can lead to non-specific signals. Try different lot of antibodies.
- Dry Plate - Do not allow the plates to completely dry during any step of the procedure. Immediately add new solutions after all washes.
- Detection Reagent concentration too low → Titrate detection reagent to higher concentration in order to increase saturation and signal.

2. High Background Signal or Non-Specific Ubiquitination

Possible Causes & Solutions:

- Insufficient washing between steps → Inadequate washing can leave non-specific proteins on the plate, leading to high background. Ensure proper plate washing using the recommended buffer and multiple wash steps.
- Excess Cell Lysate → Optimize lysate concentrations to obtain best signal to background
- Primary antibody concentration too high → Titrate the antibody to reduce non-specific binding.
- Non-specific ubiquitination of other proteins → Run a negative control (no cell lysate) to confirm specificity.
- Contaminated reagents or buffers → Use fresh reagents and ensure proper storage.
- If using a plate washer to wash the plates, use freshly made wash buffer.
- Bubble Formation - Avoid creating bubbles. Remove bubble before taking the readings as bubbles can interfere with the readings.
- Detection Reagent Concentration too high → Titrate detection reagent down in order to lessen the saturation allowing for a bigger difference between samples

3. Unexpected Ubiquitination Patterns

Possible Causes & Solutions:

- Cell lysate instability → Avoid multiple freeze-thaws to prevent unnecessary proteolysis.

K63 Ubiquitin Linkage ELISA Kit (Chain Selective)

4. Inconsistent Results Between Experiments

Possible Causes & Solutions:

- Variation in cell lysate or reagent concentrations → Always prepare master mixes and use consistent cell lysate samples.
- Temperature fluctuations → Maintain a stable incubation temperature throughout the assay.
- Improper washing of plates → Inconsistent washing can cause variability in signal intensity. Ensure uniform, thorough washing between steps.
Poor pipetting accuracy → Use calibrated pipettes and low-retention tips to ensure precise reagent additions. Use a multichannel pipette for quick and uniform addition of solution to wells.
- Cross-contamination between wells → Be cautious when adding reagents to prevent splashing or cross-contamination between wells.
- Incomplete Reactions → Ensure that the reagents and samples are added to the bottom of the well. Complete mixing of solutions in the well is critical.

5. Poor Reproducibility Across Experiments or Between Users

Possible Causes & Solutions:

- Inconsistent incubation times → Use a timer to ensure precise reaction timing between experiments.
- Variation in reagent lot numbers → If switching reagent lots, validate the new batch by running side-by-side comparisons.
- Improper mixing of reagents → Always mix reagents gently but thoroughly to avoid gradients in concentration within the wells.
- User variability in pipetting technique → Train all users on consistent pipetting methods and consider using multi-channel pipettes to minimize variability.

Final Tips for Experimental Success

- Run Controls – Always include positive and negative controls to validate results.
 - **Use High-Quality Antibodies** – Poor antibodies can produce misleading data. Choose ELISA-validated antibodies with strong supporting data.
 - Optimize Titrations – Proper titration of cell lysate, primary antibody, and detection reagent is essential for high-quality results.
 - Check Reagent Stability – Ensure cell lysate samples are stored correctly to prevent degradation.
 - Repeat Key Experiments – If results are inconsistent, repeat experiments with fresh reagents and carefully controlled conditions.
 - Wash the Plate Properly – Inadequate washing between steps can lead to high background signals and retention of non-specific proteins. Be sure to follow recommended washing steps, including multiple wash cycles, to ensure cleaner results.
-

K63 Ubiquitin Linkage ELISA Kit (Chain Selective)

LIMITED USE LABEL LICENSE

The K63-Chain Selective Assay Plate and associated technologies are licensed from LifeSensors, Inc. For information on obtaining a license for commercial purposes, contact Director of Business Development, info@lifesensors.com, LifeSensors, Inc., Malvern PA 19355, Phone: 610.644.8845 or www.lifesensors.com.

The product and/or its derivatives are being furnished to the purchaser under the following conditions: By purchasing the product, the purchaser agrees to comply with the terms of this Limited Use Label License. The product is provided to purchaser for research use only. The purchaser of this product may not transfer or otherwise sell this product or its derivatives to a third party and no rights are given to the purchaser to use the product or its derivatives for commercial purposes as defined below.

Commercial purposes mean any activity for which a party receives consideration and may include, but is not limited to,

1. use of the product or its derivatives in manufacturing,
2. use of the product or its derivatives to provide a service, information, or data,
3. use of the product or its derivatives for diagnostic purposes,
4. transfer or sale of protein or chemical component made with the product or derivatives of the product,
5. re-sale of the product or its derivatives, whether or not such product or its derivatives are resold for use in research.

Product use must follow compliance with all laws and regulations, including but not limited to current EPA, FDA, USDA, and NIH guidelines. THE MATERIALS WILL NOT BE USED IN HUMANS.

Purchaser acknowledges that the product is experimental in nature and provided WITHOUT WARRANTY OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE OR ANY OTHER WARRANTY, EXPRESS OR IMPLIED. LIFESENSORS, INC. MAKES NO REPRESENTATION OR WARRANTY THAT THE USE OF THE PRODUCT WILL NOT INFRINGE ANY PATENT OR OTHER PROPRIETARY RIGHTS.

In no event shall LifeSensors, Inc. be liable for any use of the product by the purchaser. Purchaser agrees to defend, indemnify, and hold harmless LifeSensors, Inc. its officers, employees, and agents from any loss, claim, injury, damage, expense, or liability (including attorney's fees), of whatsoever kind or nature, which may arise from or in connection with this Agreement, including but not limited to purchaser use, handling or storage of the product.

K63 Ubiquitin Linkage ELISA Kit (Chain Selective)

all products are for research use only • not intended for human or animal diagnostic or therapeutic uses
LifeSensors, Inc., 271 Great Valley Parkway, Malvern PA 19355 • (p) 610.644.8845 (f) 610.644.8616
techsupport@lifesensors.com • www.lifesensors.com • sales@lifesensors.com
Copyright © 2010 LifeSensors, Inc. All Rights Reserved