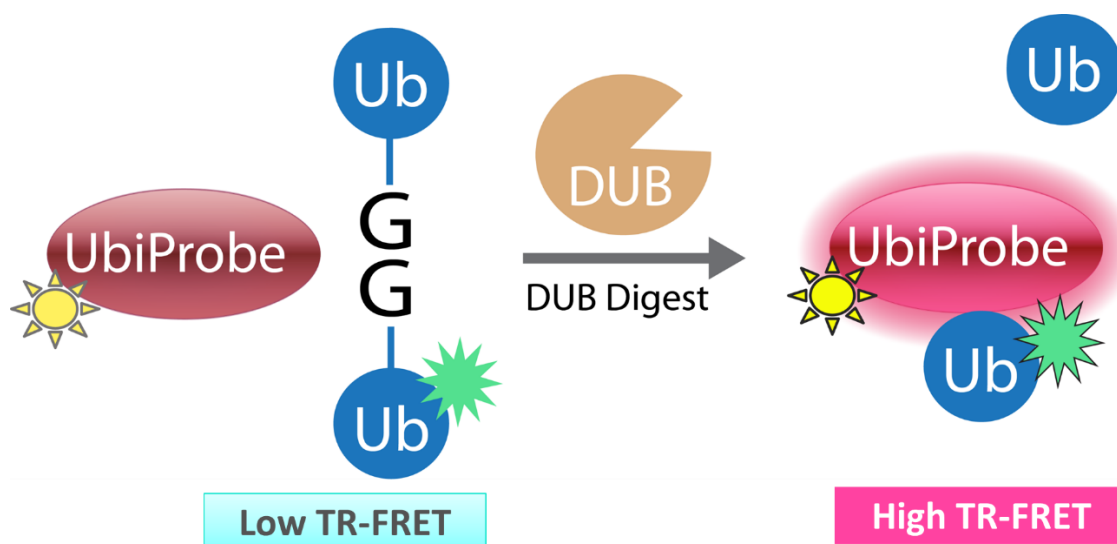


UbiProbe: K63 Di-Ub DUB Assay Kit (Catalog # [DU501](#))

MANUAL

UbiProbe: K63 Di-Ub DUB Assay Kit

Catalog Number: [DU501](#)



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

UbiProbe: K63 Di-Ub DUB Assay Kit (Catalog # [DU501](#))

UbiProbe: K63 Di-Ub DUB Assay Kit

Cat. # DU501

BACKGROUND

The UbiProbe assay consists of a designed Di-Ubiquitin (Di-Ub) fluorophore labeled as a reporter substrate, along with a ubiquitin binding reagent called "UbiProbe" serving as a detection probe. This arrangement generates a TR-FRET signal indicative of Deubiquitinase (DUB) enzyme activity. In the absence of DUB or in presence of DUB inhibitors, the uncleaved Di-Ub reporter substrate is considered inactive, leading to low TR-FRET signal. Upon cleavage of the Di-Ub-reporter substrate by the DUB, the cleaved ubiquitin interacts with UbiProbe, producing a TR-FRET signal. This coupled assay utilizes the signal generated via the cleavage of the Di-Ub substrate as a quantitative measure of isopeptidase activity in a homogeneous high-throughput format. Different Di-Ub reporter substrates (available with M1, K6, K11, K33, K48, and K63 linkages) allow for demonstration of selectivity and evaluation chain-specific DUBs.

SUGGESTED USES:

1. Real-time monitoring of DUB activity against DiUbiquitin substrates.
2. Profiling and validation of DUB inhibitors or activators.
3. High throughput screening (HTS) to identify DUB inhibitors.
4. Selectivity studies and DUBTACs screening.
5. Mechanism of action studies of DUB inhibitors or activators.

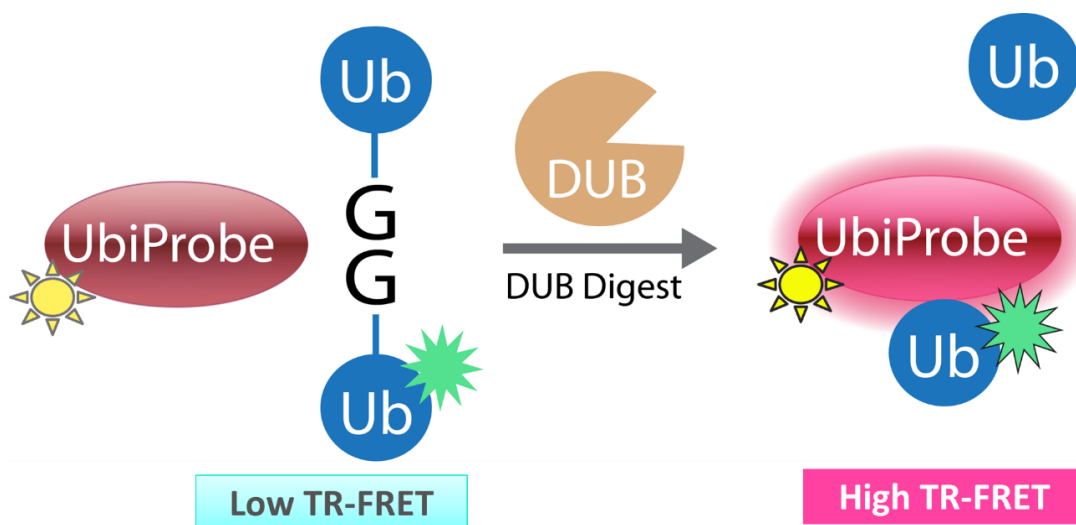


Figure 1. Schematic of UbiProbe DUB Assay. Cleavage of DiUb substrate by DUB results in high TR-FRET signal.

UbiProbe: K63 Di-Ub DUB Assay Kit (Catalog # [DU501](#))

COMPONENTS

Store all materials at -80°C, avoid cycles of freezing and thawing. All components are stable for at least 2 months.

1. **10X Assay Buffer**
Size: 1.0 mL (10X)
Note: Add β -mercaptoethanol fresh to a final concentration of 1mM in 1X assay buffer.
2. **UbiProbe**
Size: 1x 18 μ l (125X, 5 μ M)
3. **K63-DiUb^{Alexa647}**
Size: 1x 18 μ l (125X, 5 μ M)
4. **MonUb^{Alexa647}**
Size: 1x 5 μ l (250X, 5 μ M)
5. **Positive Control DUB, USP2core**
Size: 1x 5 μ l (400X, 20 μ M)

ADDITIONAL ITEMS REQUIRED BUT NOT INCLUDED IN THE KIT

1. Deubiquitinase of your choice. [DUBs](#) can also be purchased separately from [Lifesensors Inc.](#)
2. Low-speed centrifuge with a microplate adaptor.
3. Multichannel pipettes suitable for dispensing 0.5 μ l and 5 μ l volumes.
4. 384 shallow-well black assay plate such as NUNC Catalog # 267461.
5. β -Mercaptoethanol (BME).
6. Streptavidin Europium Chelate W-1024 (Sa-Eu) purchased from Columbia Biosciences (Cat# D17-2212-50).
7. TR-FRET capable plate reader with UV (320 nm +/- 20 nm), 620 nm, and 665 nm optical filters.
8. 96 or 384 well polypropylene plate for preparing compound dilutions.
9. Plate seals (optional).

IMPORTANT NOTES

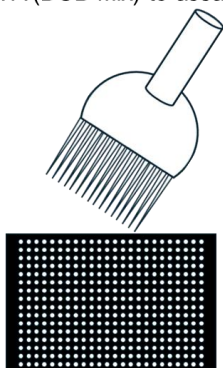
1. **Do not add 10X assay buffer directly to DUB, UbiProbe, Di-Ub Substrate or Streptavidin Europium.**
2. **After thawing the reagents, avoid making small volume (for example 2 μ l) aliquots of UbiProbe and Di-Ubiquitins for freezing at -80°C. Components have been validated for activity after multiple freeze-thaw cycles.**

UbiProbe: K63 Di-Ub DUB Assay Kit (Catalog # [DU501](#))

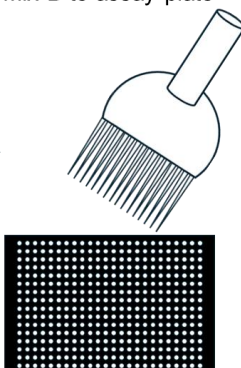
3. *Streptavidin Europium should be stored according to manufacturer recommendations, typically at 2 °C - 8 °C.*

ASSAY SUMMARY

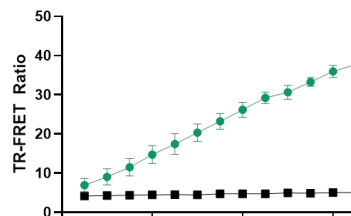
Step 1: Dispense 5 µl per well Mix A (DUB Mix) to assay plate



Step 2: Dispense 5 µl per well Mix B to assay plate



Step 3: Read Assay Plate with TR-FRET capable plate reader.



ASSAY SETUP

– See Table 1 for example concentrations and volumes.

1. **Prepare 1x Assay buffer.** Dilute 10x assay buffer in water and add β-Mercaptoethanol (BME) to a final concentration of 1.0 mM. For example, dilute 1 mL of 10x assay buffer in 9 mL water and add 0.7 µl BME (14.4 M). If testing compounds, dispense 0.5 µl of dimethylsulfoxide (DMSO) or compounds at a 21x concentration to the assay plate.

Note. A reducing agent must be added to the assay buffer for optimal DUB activity. DTT (5 to 10 mM) can be used as an alternative to β-mercaptoethanol.

2. **Prepare Mix A:** Prepare Mix A by diluting your deubiquitinase (DUB) at a 2x screening concentration in 1x assay buffer. For example, to set up a USP2Core assay with a final concentration of 5 nM, prepare a 2x solution by adding 10 nM USP2Core in the required volume of 1x assay buffer.
3. Dispense 5 µl of Mix A per well to the assay plate. Include blank control wells and positive control wells to ensure that the assay is working properly.
 - a. For a blank control (-DUB), dispense 5 µl of assay buffer into corresponding wells.
 - b. For a positive control, prepare a 40 nM solution of MonoUb^{Alexa647}. Dispense 5 µl/well. The MonoUb^{Alexa647} will bind UbiProbe and generate a high TR-FRET signal.
4. Briefly centrifuge assay plate for ~30 seconds at ~2000 rpm. If testing compounds, the DUB may be preincubated with the compounds for 30min at room temperature. Seal or cover the assay plate until the next step.
5. **Prepare Mix B:** Prepare Mix B by diluting UbiProbe, K63-DiUb^{Alexa647}, and Streptavidin Europium in assay buffer at 2x concentrations. For example, prepare a 2x mix containing 40 nM UbiProbe, 40 nM K48-DiUb^{Alexa647} substrate and 2 nM Streptavidin Europium in 1x assay buffer. Dispense 5 µl per well to the assay plate.

UbiProbe: K63 Di-Ub DUB Assay Kit (Catalog # [DU501](#))

6. Briefly centrifuge assay plate for ~30 seconds at ~2000 rpm.
7. Read the assay plate in a plate reader capable of measuring TR-FRET signal. See table 2 for suggested plate reader parameters. For DUBs that are highly active or DUBs at high concentration (for example, >100 nM), assay kinetics may be very rapid. It is recommended to titrate the DUB concentration to identify the optimal concentration that exhibits linear assay kinetics.

EXAMPLE CONCENTRATIONS AND VOLUMES

Table 1. Example Volume calculations for a single full 384-well plate

Mix A (2.25 ml)	2x Concentration	Volume	Final concentration
Assay Buffer	1x	2.25 ml	1x
USP2c (20µM)	10 nM	1.13 µl	5 nM
Positive TR-FRET Control (125 µl)			
Assay Buffer	1x	125 µl	1x
Mono-Ub ^{Alexa647} (5 µM)	40 nM	1 µl	20 nM

Mix B (2.25 ml)	2x Concentration	Volume	Final concentration
Assay Buffer	1x	2.25 ml	1x
UbiProbe (5 µM)	40 nM	18 µl	20 nM
K63-DiUbiquitin-Ub ^{Alexa647} (5 µM)	40 nM	18 µl	20 nM
Streptavidin Europium (1.85 µM)	2 nM	2.43 µl	1 nM

PLATE READER SETTINGS

Table 2. Suggested plate reader parameters for Time-Resolved FRET detection

Excitation	Wavelength	Integration Time Start	Integration Time Stop
Excitation	320 nm, ± 10 nm		
Channel 1 Emission	620 nm, ± 10 nm	60 µs	400 µs
Channel 2 Emission	665 nm, ± 10 nm	60 µs	400 µs

UbiProbe: K63 Di-Ub DUB Assay Kit (Catalog # [DU501](#))

REPRESENTATIVE DATA

1. Validation of USP2core as a pan-linkage specific DUB using UbiProbe kits

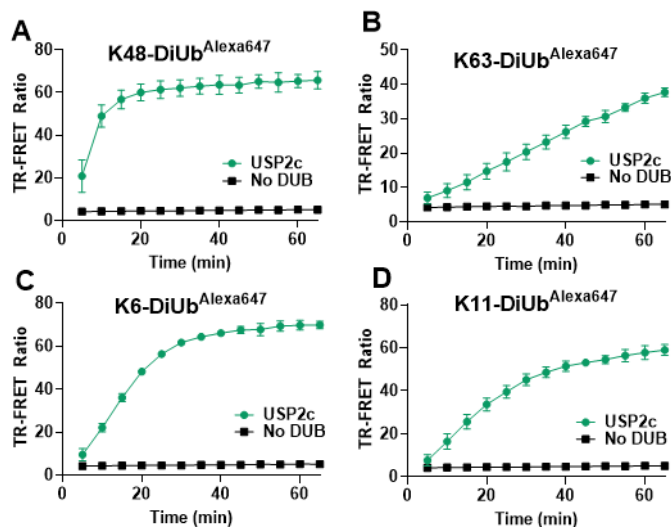


Figure 2. Robust USP2 activity obtained using UbiProbe assay with various DiUb substrates. [USP2Core](#) (5 nM) was incubated with [K48-DiUb](#) (A), [K63-DiUb](#) (B), [K6-DiUb](#) (C) and [K11-DiUb](#) (D) and TR-FRET was measured for 65 minutes at 5-min intervals. DUB activity with each of the Di-ubiquitin probes is represented as TR-FRET ratio and compared to negative control (-DUB). The TR-FRET signal was acquired using BMG Labtech ClarioStar. The error bars represent standard deviation, n=4.

UbiProbe: K63 Di-Ub DUB Assay Kit (Catalog # [DU501](#))

REFERENCES

1. Hershko, A. and A. Ciechanover, The ubiquitin-proteasome pathway. *Annu Rev Biochem*, 1998. **67**: p. 425-479.
2. Schauer NJ, Magin RS, Liu X, Doherty LM, Buhrlage SJ. Advances in Discovering Deubiquitinating Enzyme (DUB) Inhibitors. *J Med Chem*. 2020 Mar 26;63(6):2731-2750. doi: 10.1021/acs.jmedchem.9b01138. Epub 2019 Nov 20. PMID: 31682427.
3. Mevissen TET, Komander D. Mechanisms of Deubiquitinase Specificity and Regulation. *Annu Rev Biochem*. 2017 Jun 20;86:159-192. doi: 10.1146/annurev-biochem-061516-044916. Epub 2017 May 12. PMID: 28498721.

UbiProbe: K63 Di-Ub DUB Assay Kit (Catalog # [DU501](#))

LIMITED USE LABEL LICENSE

The UbiProbe: K63 Di-Ub DUB Assay Kit 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.