

Ubiquitin Conjugating Enzymes and Conjugation Assays: Deciphering the Complexity of Ubiquitylation

We offer a large selection of enzymes to study the multi-step process of ubiquitylation: E1-E2-E3.

It is as easy as 1-2-3!

Create your own assay

OR

Use our preconfigured ubiquitin conjugation assays to measure the activity of ubiquitin conjugating enzymes

- Microtiter plate-based
- High Throughput Screening
- A wide range of E2s and E3s

APPLICATIONS:

Quantitation of E2 and/or E3 Ligase activity

Demonstration of novel E2 or E3 Activity

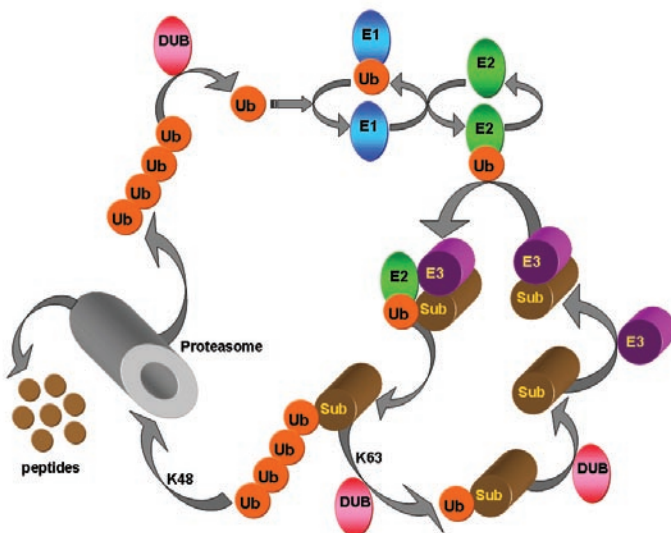
Identification of novel substrates for E3 of Interest

Confirmation of Putative E3 substrates

High throughput screening (HTS) for antagonists or agonists of ligases

Monitoring enzyme activity during purification

E1-E2-E3 conjugation machinery



Ref: Marblestone et al, 2010 (in press)

Contact Us
To Learn More:
info@lifesensors.com

To order, please call
 610-644-8845 or visit
www.lifesensors.com

Select a single enzyme for your experimental design or create a customized assay

E1, E2, E3 – see our website to make a selection

E2 customized profiling kit or E2 selectivity panel –select from over 25 ligases and save \$\$!

NEW

E₃LITE Custom Ubiquitin Ligase Kit

NEW

E3 Substrate Identification Service and Kit for Microarrays: **DIY or our team will do it for you!**

Novel, easy to customize tools to study the complicated world of ligases:

Create an assay that meets your expectations

E₃LITE Custom Ubiquitin Ligase Kit

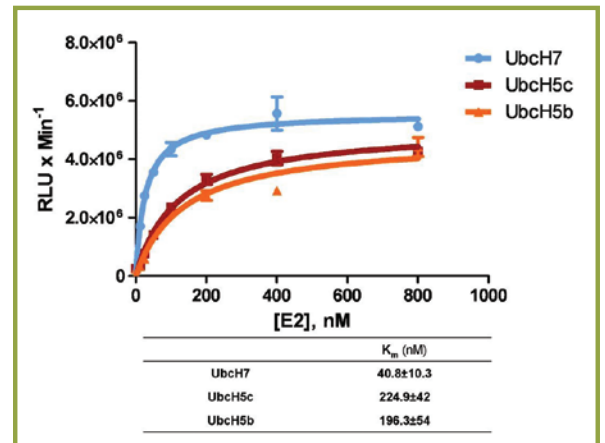
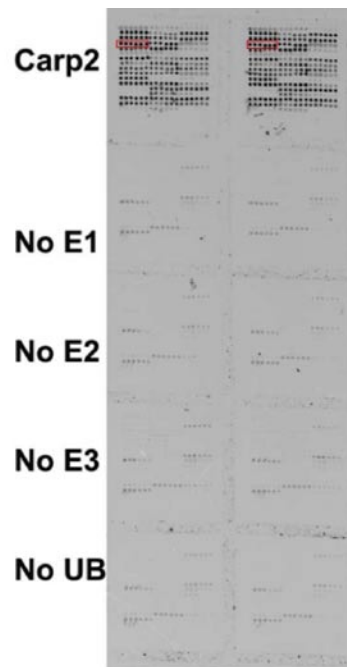
- Quantitative measure of E3 ligase activity
- Choice of E2
- Native, unlabeled ubiquitin

E3 Substrate Identification Kit for Microarrays

- Identification of novel substrates for E3 enzymes
- Novel, sensitive poly-ubiquitin detection reagent with nanomolar affinity
- Over 20 E2 conjugation enzymes representing members of each enzyme class

Generate reliable data

Measuring Km values for E2 conjugating enzymes with the E3 ligase, E6AP



E3 Substrate Identification Kit for microarrays example

