Mix-and-Read Assays For Regulated Ubiquitination & Degradation of Proteins Using PathHunter® Technology

Daniel Bassoni, William Raab, Mong Saetern and Tom Wehrman, Ph.D. DiscoveRx Corporation, Fremont, CA 94538, USA









Both β-catenin and NRF2 are regulated by continuous ubiquitination and degradation in the absence of activating stimuli by the β-TRCP/SCF and KEAP1-cul3 complexes respectively. In the presence of signal the proteins are stabilized and translocate to the nucleus. To monitor this process, cells were engineered to express the target fused to the ePL tag, with the second complementing partner (EA) expressed in the nucleus. Upon stabilization and translocation of the ePL tagged protein, complementation ensues and signal is detected using the PathHunter detection reagents. The top panel shows U2OS cells that are expressing β-catenin fused to ePL are stimulated by a titration of Wnt3A to activate endogenous Frizzled receptors, and subsequently, β-catenin. Cells were incubated with ligand for 6 hours and signal was detected using Pathhunter detection reagents. In the bottom panel, cells expressing NRF2-ePL, KEAP1, and EA (in the nucleus) were incubated for 6 hours in the presence of antioxidant compounds. Signal was also detected using the PathHunter detection reagents.

Summary of the ePL tag advantages for ubiquitination and regulated proteolysis

- Minimally invasive tag = natural protein life cycle -
- Sensitivity to low pM levels levels of the fusion proteins to be detected.
- A homogenous, whole cell, single addition assay
- Ready to go Biosensor cells
- Highly flexible -



PathHunter[®] Pathway Assays Ready-to-go, biosensor cell lines with a complimentary detection system offers you convenience and saves months of assay development time.

Custom Projects Utilize DiscoveRx's proprietary EFC technology to build your own functional cell-based kinase assays. Talk to our experts about your kinase targets, contact CAD@discoverx.com.

Custom Screening & Profiling Services Send your compound for family-wide profiling in select targets or for selectivity studies and unique information. For more information, please email profiling@discoverx.com.

© 2013 DiscoveRx Corporation. All Rights Reserved.

EFC APPLICATION TO OTHER UBIQUITIN REGULATED PATHWAYS

SUMMARY

- This maximizes the assay window by not imparting stability or instability from the detection tag.

- Each ePL fusion creates an enzyme, upon addition of the InCELL detection reagent, allowing low

- Highly miniaturizable and readily adaptable, from bench-top to HTS.

- Assay ready kits, clonal cell lines, and custom assay options for your target of interest.

- N and C-terminal fusions perform comparably, the ePL can be used by itself for degradation assays or in conjunction with EA for translocation assays.

TECHNOLOGY ACCESS

DRX POSTER PH-TECH UBIOUITIN 0213

