

Cellular-HTS Assays for the Ubiquitin-Proteasome system in NF κ B Signaling

Matthew Robers, Sara Hereley, Robert Horton, and Thomas Machleidt
Invitrogen Corporation • 1600 Faraday Avenue • Carlsbad, California 92008 • USA

Introduction

Activation of the NF κ B pathway induces a signal transduction cascade that results in the phosphorylation, ubiquitination, and proteasomal degradation of I κ B α . Upon I κ B α degradation, liberated NF κ B translocates to the nucleus to activate target gene expression (Fig 1). We have developed a set of target-specific HTS-compatible assays capable of interrogating the ubiquitin-proteasome system. Using a clonal cell line expressing GFP-I κ B α , we have developed a TR-FRET assay capable of measuring ubiquitination of endogenously expressed I κ B α . We have concurrently developed a living-cell β -lactamase (*bla*) reporter assay for degradation of Bla-I κ B α fusion proteins. These technologies provide a powerful means to interrogate the intermediate steps in NF κ B signaling, without compromising the endogenous physiological complexity of this signaling pathway. Cellular HTS assays that interrogate these processes will provide a unique integrated approach to dissecting the ubiquitin-proteasome system in the context of NF κ B signaling.

Figure 1 – NF κ B Pathway

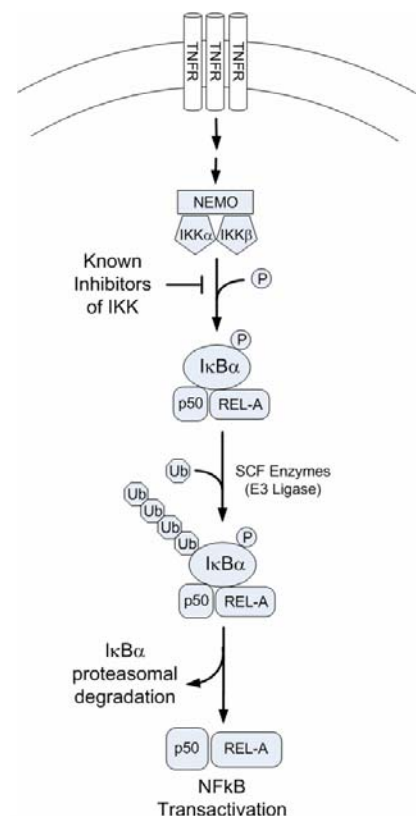


Figure 2 – Inducible degradation of GFP-I κ B α in HEK-293 cells by Western analysis



Figure 2 A clonal GFP-I κ B α -expressing cell line was generated in order to probe the post-translational modifications of I κ B α . We validated this construct for the inducible depletion of GFP-I κ B α (approximately 65 kDa) by treating the cells with TNF α , generating lysates from each sample, and performing Western Analysis using anti-GFP antibodies.

Figure 3 – Assay schematic for ubiquitination of GFP-I κ B α expressed in HEK-293 cells

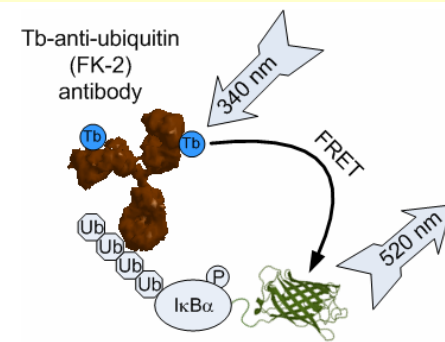


Figure 3 In cell lysates, Tb-anti-ubiquitin antibodies bind ubiquitinated GFP-I κ B α , allowing the Tb and GFP fluorophores to come in close proximity for energy transfer to occur.

Figure 4 – Workflow for cellular GFP-I κ B α ubiquitination assay

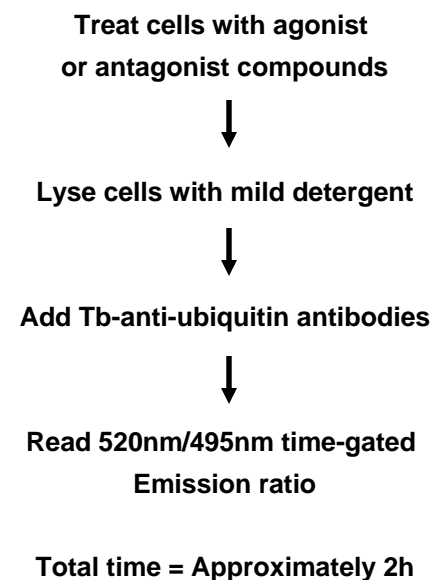
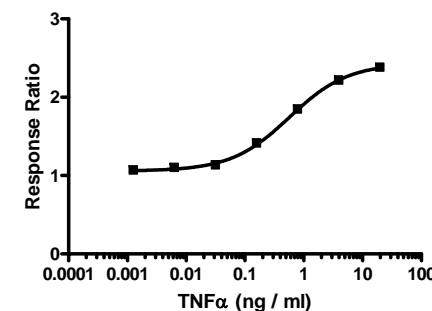


Figure 5 – Lanthascreen™ TR-FRET assays for ubiquitination of GFP-I κ B α in HEK-293 cells

A TNF α -induced ubiquitination of GFP-I κ B α



B Inhibition of ubiquitination of GFP-I κ B α

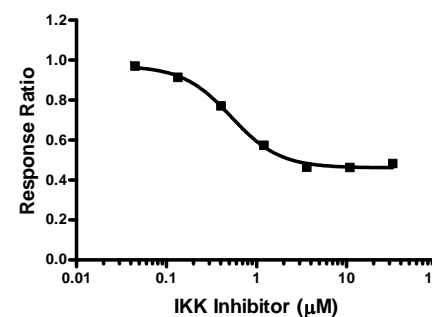
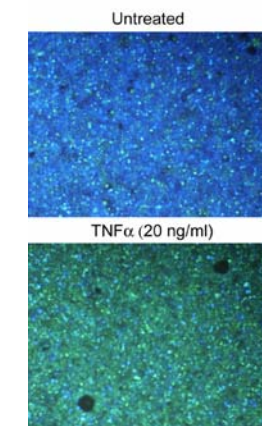


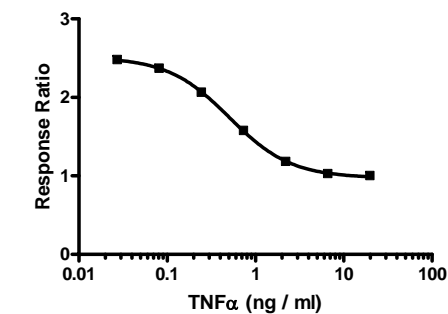
Figure 5 a. In cell lysates, a dose-dependent rise in TR-FRET signal is seen, concomitant with an increase in ubiquitinated GFP-I κ B α . Briefly, in 96 well format this cell line was treated with serial dilutions of TNF α , lysates were generated from each sample, and the lysate was probed with 10nM Tb-anti-ubiquitin antibody solution. A time-gated fluorescence emission ratio was then generated for each sample using the 520 nm (GFP) emission signal referenced against the 495 nm (Tb) emission signal. Response ratios were generated by dividing each emission ratio value by the unstimulated value (zero agonist or antagonist). Measurements were taken using a Tecan Ultra Fluorescence plate reader. **b.** Inhibition of TNF α -induced ubiquitination using I κ B-kinase inhibitor IV. This cell line was pretreated with serial dilutions of inhibitor compound prior to stimulation with 1ng/mL TNF α .

Figure 6 – β -Lactamase assays for Bla-I κ B α degradation in living HEK-293 cells

A Image Analysis of Bla-I κ B α degradation



B TNF α -induced degradation of Bla-I κ B α



C Inhibition of degradation of Bla-I κ B α

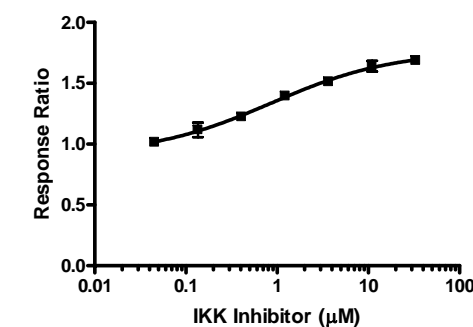


Figure 6 a. Color FRET images of Bla-I κ B α degradation using the cell-permeable substrate for Bla, CCF4-AM. When intracellular Bla levels are above a threshold, the green-emitting CCF4-AM substrate is cleaved to generate a blue product. Thus, untreated (blue) cells become green in response to TNF α . **b.** In living HEK-293 cells, a dose-dependent depletion of Bla activity is seen, concomitant with the proteolytic processing of Bla-I κ B α . FRET emission ratios were generated using the 460 nm value (blue) divided by the 530 nm value (green). Response ratios were generated by dividing each blue/green ratio value by the unstimulated value (zero agonist or antagonist). **c.** Inhibition of TNF α -induced degradation using I κ B-kinase inhibitor IV. This cell line was pretreated with serial dilutions of inhibitor compound prior to stimulation with 1ng/mL TNF α . Data was generated using a Tecan Safire II fluorescent plate reader.