

EZClick™ Global Protein Synthesis Assay Kit (FACS/Microscopy), Green Fluorescence rev 06/20

(Catalog # K459-100; 100 assays; Store at -20°C)

I. Introduction:

Cells generate a complete set of proteins during division. Protein synthesis is a tightly regulated process. Many critical controls in gene expression occur at the translational level to ensure that production of specific cellular proteins is quickly turned on/off under specific conditions (heat shock, starvation, etc.). Protein synthesis is essential for cell growth, proliferation, signaling, differentiation or death. Therefore, the identity and amount of the synthesized proteins are critical parameters in determining the physiological state of the cell. Methods enabling the detection and characterization of nascent proteins, or changes in spatial and temporal protein expression/degradation patterns during disease, drug treatments or environmental changes are important tools in assessment of cytotoxicity. **BioVision's EZClick™ Global Protein Synthesis Assay Kit** utilizes a novel and robust chemical method based on an alkyne containing and cell-permeable analog of puromycin, O-Propargyl-puromycin (OP-puro). Once inside the cell, OP-puro stops translation by forming covalent conjugates with nascent polypeptide chains. Truncated polypeptides are rapidly turned over by the proteasome and can be detected based on a click reaction with the fluorescent azide. Unlike methionine analogs, OP-puro does not require methionine-free conditions and can be used to label nascent proteins directly in the cell culture. Our kit provides sufficient materials for 100 assays to detect nascent proteins synthesized under various physiological conditions, and Cycloheximide, an inhibitor of protein synthesis that serves as an experimental control.

II. Applications:

- Detection of nascent protein biosynthesis
- Detection of protein expression or degradation patterns in presence of cytotoxic agents
- Screening for genotoxic compounds and effectors of protein synthesis

III. Sample Type:

- Suspension or adherent cell cultures

IV. Kit Contents:

Components	K459-100	Cap Code	Part Number
EZClick™ Wash Buffer (10X)	25 ml	NM	K459-100-1
Fixative Solution	10 ml	WM	K459-100-2
Permeabilization Buffer (10X)	25 ml	NM/Blue	K459-100-3
EZClick™ Protein Label (400X)	25 µl	Clear	K459-100-4
Copper Reagent (100X)	100 µl	Blue	K459-100-5
EZClick™ Fluorescent Azide (100X)	100 µl	Green/Amber Vial	K459-100-6
Reducing Agent (20X)	500 µl	Yellow	K459-100-7
EZClick™ Total DNA Stain (1000X)	20 µl	Blue/Amber Vial	K459-100-8
Cycloheximide (100X)	10 µl	Green	K459-100-9

V. User Supplied Reagents and Equipment:

- Tissue culture vessels and appropriate culturing media
- Phosphate Buffered Saline (PBS, pH 7.4)
- Sterile 0.1% Gelatin Solution (optional, only required for suspension cells)
- Flow cytometer equipped with laser capable of excitation at 488 nm wavelength (FL-1)
- Fluorescence microscope capable of excitation and emission at 440/490 and UV filter

VI. Storage Conditions and Reagent Preparation:

Upon arrival, store the entire kit at -20°C protected from light. Briefly centrifuge small vials prior to opening. Read the entire protocol before performing the assay.

- **10X Wash Buffer and 10X Permeabilization Buffers:** Thaw at 37°C to dissolve completely. Dilute the 10X stocks 1:10 in sterile water, mix well. Store at 4 °C.
- **Fixative Solution:** Ready to use. After opening, store at -20°C, protected from light.
- **Remaining components:** Store at -20°C protected from light. While in use, keep on ice and minimize light exposure.

VII. Assay Protocol:

Notes:

This assay was developed with HeLa (adherent) and Jurkat (suspension) cells and can be modified for any cell line. The protocol below refers to a 96-well tissue culture plate; adjust volumes accordingly for other plate formats. The assay volume is 100 µl. Growth conditions, cell number per well and other factors may affect the incorporation rate of the Protein Label; therefore optimize the assay for your cell type. We suggest an initial test of several EZClick™ Protein Label concentrations to find best conditions for your experimental design. Avoid stressing the cells by washes or temperature changes prior to incubation with EZClick™ Protein Label. All steps should be carried out at room temperature (RT) unless otherwise specified; equilibrated all buffers to RT prior to the experiment

1. Labeling of control and experimental cells; method with **drug pre-incubation**:

- Obtain cell suspension of desired density and seed directly into tissue culture vessels, or on coverslips for high resolution microscopy. **To immobilize suspension cells for microscopy:** add 100 µl of 0.1% gelatin solution directly into the wells. Tilt the plate to cover the entire well surface and place it in a tissue culture hood for 1 hour. Gently remove the gelatin solution and seed your cells. Allow the cells to recover overnight before the treatment. Next day, treat the cells with appropriate effectors according to your protocol; *do not add treatment to the positive and negative control cells.* **Negative Control** (cells not exposed to the Protein Label or treatment), **Positive Control** (cells incubated with 1X Protein Label only).
- To use Cycloheximide as an inhibitor of protein synthesis, dilute it 1:100 in the culture medium and incubate the cells for 30 minutes at a 37°C. Remove the drug-containing media. **For immobilized suspension cells:** Centrifuge the plate at 300 x g (or the lowest centrifuge setting) for 5 minutes to deposit the cells onto the surface. Tilt the plate and **gently** remove the media with a pipette tip. It is important to avoid excessive centrifugation speeds, which can damage the cells. *Make note of the place that is used, and perform subsequent aspirations from the same place.*

FOR RESEARCH USE ONLY!

- c. Replace the media with fresh aliquots containing EZClick™ Protein Label (400X) diluted to 1X final concentration and tested compound or Cycloheximide. Add Protein Label to the positive control cells. Incubate the cells for additional 0.5 - 2 hours in a 37°C incubator, or for the period of time required by your experimental protocol. **Note:** for drug and EZClick™ Protein Label **co-incubation**, dilute EZClick™ Protein Label directly into the drug or Cycloheximide treated cells, do not change the media.
- d. Terminate the experiment by removal of the culture medium and rinse the cells once with 100 µl of PBS, discard the supernatant. Proceed to the Fixation and Permeabilization.

2. Fixation and Permeabilization:

- a. **For adherent cells:** Add 100 µl of Fixative Solution to each well and incubate the cells for 15 min at RT protected from light. Remove the fixative and wash the cells once with 200 µl of 1X Wash Buffer, remove the wash. Add 100 µl of 1X Permeabilization Buffer and incubate the cells for 10 min at RT. Remove the Permeabilization Buffer.
- b. **For suspension cells:** Re-suspend the cells in 100 µl of Fixative Solution and incubate for 15 min at RT protected from light. Centrifuge cells at 900 x g for 5 minutes and remove the fixative Solution. Wash the cells once with 200 µl of 1X Wash Buffer. Centrifuge cells at 900 x g for 5 minutes and discard the supernatant and re-suspend the cells in 100 µl of 1X Permeabilization Buffer. Incubate the cells for 10 min at RT. Remove the Permeabilization Buffer.

3. EZClick™ Protein reaction and total DNA staining:

- a. **Reaction cocktail:** Prepare 1X EZClick™ reaction cocktail according to the table below. Volumes should be multiplied by number of samples and reagents added in the exact order. Use the reaction cocktail within 15 minutes of preparation. *Cells should be protected from light during, and following the EZClick™ reaction and DNA staining.*

	Amount per Reaction
PBS	93 µl
Copper Reagent (100X)	1 µl
EZClick™ Fluorescent Azide (100X)	1 µl
Reducing Agent (20X)	5 µl

- b. **EZClick™ Protein Reaction:** Add 100 µl of 1X EZClick™ Reaction cocktail to each sample (Negative Control Cells: Add 100 µl PBS only) and incubate the cells for 30 min at RT protected from light. Centrifuge cells at 900 x g for 5 minutes and remove the reaction cocktail and wash cells in 200 µl of Wash Buffer. Centrifuge cells at 900 x g for 5 minutes and remove the wash buffer. Do 2 times Wash Step and suspend the cells in 100 µl of PBS. Proceed to DNA staining. If no DNA staining is desired, proceed to Microscopic or FACS analysis. **DNA staining:** Prepare 1X dilution of Total DNA Stain and add 100 µl per well. Incubate the cells for 20 minutes at RT, or refrigerate at 4 °C protected from light. Centrifuge cells at 900 x g for 5 minutes and remove the DNA stain solution; wash the cells once with 100 µl of PBS. Note: cells are compatible with all methods of slide preparation including wet mount or prepared mounting media.
4. **Fluorescence Microscope Imaging:** Analyze samples for green fluorescence generated by labeled Protein and for blue fluorescence by nuclear DNA. **FACS analysis:** Harvest the cells by preferred method and wash with 0.5 ml of ice-cold PBS. Centrifuge cells at 900 x g for 5 minutes and re-suspend the pellets with 100 µl of ice-cold PBS and analyze samples for green fluorescence generated by *de novo* synthesized protein during click reaction.

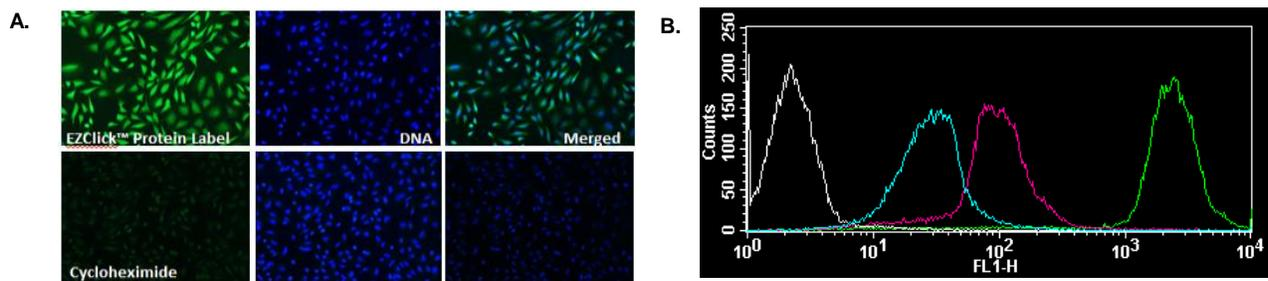


Figure: Inhibitory effect of Cycloheximide on nascent polypeptides synthesis. HeLa (10^5 cells/ml) and Jurkat (1×10^6 cells/ml) cells respectively were pre-treated with vehicle or Cycloheximide for 30 min at 37°C. Subsequently, cells were incubated for additional 30 min with fresh aliquots of media containing either EZClick™ Protein Label or EZClick™ Protein Label and Cycloheximide. Cells were then processed and analyzed by Microscopy and FACS according to the kit protocol. **(A)** Green fluorescence (upper panel) corresponds to *de novo* synthesized polypeptides whereas bottom panel shows the inhibitory effect of Cycloheximide on protein biosynthesis. Nuclear staining in both panels confirms that green signal is a result of Protein Label incorporation. **(B)** FACS analysis of negative control (white), background (Azide only, blue), positive control (Protein Label, green) and Cycloheximide-treated (pink) cell populations. Signal measured in FL-1 channel clearly shows the inhibitory effect of Cycloheximide on nascent polypeptides synthesis.

VIII. RELATED PRODUCTS:

- EZClick™ Global RNA Synthesis Assay Kit (FACS/Microscopy), Red Fluorescence (K718)
- EZClick™ Global Phospholipid Synthesis Assay Kit (FACS/Microscopy), Red Fluorescence (K717)
- EZClick™ EdU Cell Proliferation/DNA Synthesis Kit (FACS/Microscopy), Red Fluorescence (K946)

FOR RESEARCH USE ONLY! Not to be used on humans.