

Autophagy Analysis Using Object Spot Counting

Using Agilent BioTek Gen5 microplate reader and imager software to analyze the size and number of autophagosomes per nuclei



Authors

Sarah Beckman, PhD
Agilent Technologies, Inc.
Winooski, VT, USA

Nadia A. Rana, PhD
and Wini Luty
Enzo Life Sciences, Inc.
Farmingdale, NY, USA

Abstract

Autophagy is critical for the maintenance of cellular homeostasis. However, dysregulated autophagy can lead to death of healthy cells and survival of cancerous cells. This application note describes the use of CYTO-ID Autophagy detection kit in combination with automated object-based spot counting to quantitatively assess the effects of starvation and rapamycin on cellular autophagy by determining the size and number of autophagosomes per cell.

Introduction

A constant supply of nutrients is required during development to provide the energy necessary for growth, metabolism, and survival. Eukaryotic cells have evolved a variety of mechanisms to adjust their metabolic activities in response to changes in nutrient levels. Nutrient starvation, stress, or reduced availability of growth factors induces eukaryotic cells to adjust their metabolism in order to survive.¹ One of the key responses to such a stress is autophagy.

Autophagy or "self-eating" is a highly conserved process by which cells break down their intracellular components.² In a healthy cell under physiological conditions, autophagy is protective. In fact, autophagy plays a variety of important roles including maintenance of the amino acid pool during starvation, damaged protein and organelle turnover, prevention of neurodegeneration, tumor suppression, cellular differentiation, clearance of intracellular microbes, and regulation of innate and adaptive immunity.³

The first step of autophagy is formation of the phagophore, a cup-shaped double membrane. The edges of this membrane elongate and engulf portions of the cytoplasm, including intracellular material such as damaged organelles and misfolded proteins.⁴ The isolation membrane expands and its open ends fuse to form a double-membrane structure called the autophagosome. Autophagosomes then fuse with lysosomes to form autolysosomes and the contents inside the autophagosome are degraded by lysosomal hydrolases. The intracellular material is then recycled back into the cytosol.⁵

One of the most well-known inducers of autophagy is starvation. Through autophagy, amino acids and other nutrients are recycled from long-lived proteins, organelles, and other components of the cytoplasm, providing an internal reserve of nutrients. Starvation rapidly induces autophagy, in part by inactivation of the mTOR (mammalian target of rapamycin) substrate S6K.⁶ In a nutrient-rich environment, mTOR inhibitors such as rapamycin can induce autophagy.

Autophagy plays a role in both the pathogenesis and prevention of disease.⁷ This is especially true in cancer, where elimination of damaged intracellular components through autophagy suppresses tissue injury and tumor initiation. However, in an established tumor, autophagy promotes cancer progression by providing substrates for metabolism, maintaining functional mitochondria, and fostering survival.⁸

Traditional methods of autophagy analysis include electron microscopy and western blot analysis of LC3-II. Electron microscopy is limited by the necessity of specialized expertise, and open to when identifying an autophagosome structure.⁹ Furthermore, flow cytometry or western blot measurements of LC3-II do not always correlate with formation of autophagosomes and do not give per-cell numbers of autophagosomes.¹⁰ This application note describes the process of using CYTO-ID Autophagy detection dye combined with Agilent BioTek Gen5 3.03 software to analyze the effects of serum starvation and rapamycin on autophagosome number in HeLa cells. An analysis was performed with Gen5 3.03 with object spot counting capability, which allows us to determine the number of autophagosomes per cell as well as their size.

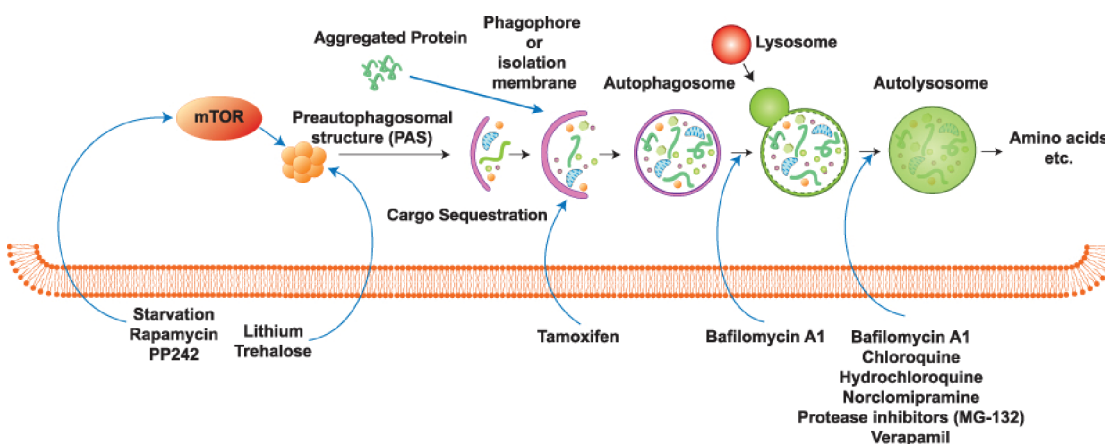


Figure 1. Schematic depiction of autophagy.

Materials and methods

Cell culture

HeLa cells were grown in Advanced Dulbecco's modified Eagle's medium (DMEM) (Gibco, Grand Island, NY) with 10% FBS (Gibco) and 1x pen/strep-glutamine (Cellgro, Manassas, VA). Cells were seeded into black-sided clear-bottom 96-well microplates from Corning (Corning, NY) at 20,000 cells per well.

Cyto-ID Autophagy detection kit

The CYTO-ID Autophagy detection kit donated by Enzo Life Sciences (Farmingdale, NY) was used to assess autophagy levels in HeLa cells. The probe is a cationic amphiphilic tracer (CAT) dye that rapidly partitions into cells. The dye is taken up by passive diffusion across the plasma membrane bilayer, and includes titratable moieties specific for selectively staining autophagic vesicles. HeLa cells were grown in normal media, which was replaced with serum-free media containing 10 μ M chloroquine for two to six hours to induce autophagy through serum starvation. Alternatively, cells were treated with 0.1 to 10 nM rapamycin with or without 10 μ M chloroquine for 18 hours to induce autophagy through mTOR inhibition. Following treatment, cells were washed 2x with 200 μ L assay buffer (1x buffer + 5% FBS). Next, the assay buffer was replaced with 100 μ L dual color detection solution (1 mL assay buffer + 1 μ L Hoechst and 2 μ L CYTO-ID green detection reagent) for 30 minutes at 37 °C in the dark. Finally, the cells were washed with 2x 200 μ L of assay buffer and this was removed and the sample was imaged in 100 μ L of assay buffer directly following the wash.

Cell imaging

Images were acquired using a 20x objective on an Agilent BioTek Lionheart FX automated microscope configured with DAPI and GFP light cubes. The DAPI light cube is configured with a 377/50 nm excitation filter and a 447/60 nm emission filter. The GFP light cube uses a 469/35 nm excitation filter and a 525/39 nm emission filter.

Image analysis

Image preprocessing was used to ensure the best possible detection of nuclei and the best separation between individual autophagosomes. Imaging preprocessing parameters are described in detail in Table 1. The GFP channel of all the images was preprocessed with a 0.5 μ m rolling ball to obtain the best separation between individual spots (Figure 2). Image preprocessing should be optimized on a per-experiment basis depending upon the size of particles being analyzed and how spread apart they are.

Table 1. Agilent BioTek Gen5 microplate reader and imager software settings. Image analysis parameters for generating a cellular mask in the DAPI channel and an object mask in the GFP channel to count autophagic vesicles.

Imaging Preprocessing	
Image Set	DAPI
Background	Dark
Rolling Bar Diameter	Auto
Image Smoothing Strength	5
Image Set	GFP
Background	Dark
Rolling Bar Diameter	0.5 μ m
Priority	Fine results
Image Smoothing Strength	0
Cellular Analysis	
Detection Channel: Primary Mask and Count	Tsf[DAPI 377,447]
Threshold	7,000
Secondary Mask	Tsf[GFP 469,525]
Measure within a Secondary Mask	Checked
Expand Primary Mask	30 μ m
Threshold	Unchecked
Count Spots	Checked
Size	0.5 to 3 μ m
Advanced Options	Count Spots Options
Rolling Ball Size	Default
Threshold	1,200

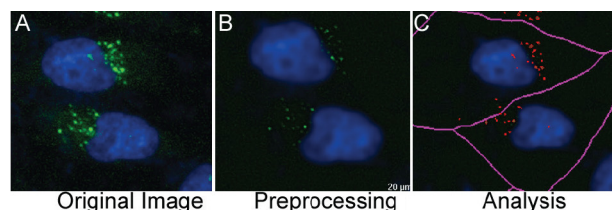


Figure 2. Autophagy spot counting workflow. (A) Original image. (B) Image after preprocessing. (C) Object masks highlighting cell area in purple and spots in red.

Object smoothing of five cycles was applied to the DAPI channel to facilitate clean masking of the nuclei. Cell counting analysis was applied to the transformed DAPI channel to highlight each individual cell. Next, object spot counting was performed on the GFP channel to determine the size and number of autophagy positive spots per nuclei according to the parameters outlined below in Table 1.

Results and discussion

HeLa cells were treated with 0.1 to 10 nM rapamycin for 18 hours to determine the effect of increasing concentration of rapamycin on the number of autophagy positive vesicles per cell. Rapamycin is an mTOR inhibitor that regulates cell growth and metabolism in response to environmental cues. Rapamycin induces autophagy due to the fact that inhibition of mTOR mimics cellular starvation by blocking signals required for cell growth and proliferation.¹ Increased CYTO-ID fluorescence indicated autophagosome formation in HeLa cells treated with rapamycin (Figure 3). There is an increase in both the size and the number of autophagic vesicles per cell as a result of increasing rapamycin concentrations (Figure 3).

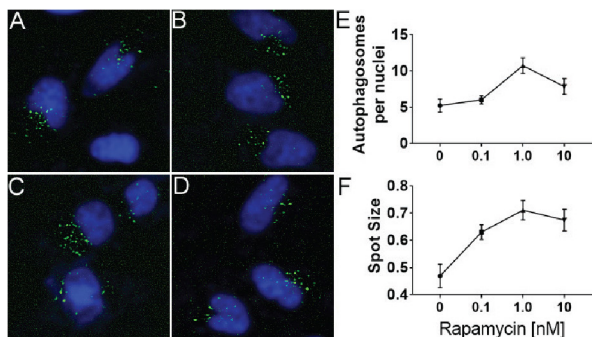


Figure 3. Autophagy-positive spot counts and size increase after treatment with rapamycin. (A) Control (B) 0.1 nM rapamycin (C) 1 nM rapamycin (D) 10 nM rapamycin (E) Autophagosomes per nuclei increase with increasing concentration of rapamycin (F) Autophagosome spot diameter increases with increasing concentrations of rapamycin.

An accumulation of autophagosomes may be indicative of either the increased generation of autophagosomes or a block in autophagosome maturation and completion of the autophagic pathway. Chloroquine is a lysosomal inhibitor that increases the pH of the lysosome, therefore preventing the activity of lysosomal acid proteases and causing autophagosomes to accumulate in the cell.¹¹ Figure 4 demonstrates an increase in the size and number of autophagy-positive spots per nuclei in response to combined rapamycin and chloroquine treatment. Notice the increase in both size and number of autophagosomes compared to rapamycin treatment alone (Figures 3 and 4).

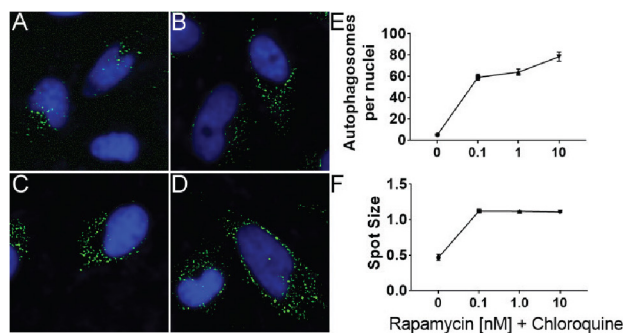


Figure 4. Autophagy-positive spot counts and size increase after treatment with rapamycin + chloroquine. (A) Control (B) 0.1 nM rapamycin + 10 μ M chloroquine (C) 1 nM rapamycin + 10 μ M chloroquine (D) 10 nM rapamycin + 10 μ M chloroquine (E) Autophagosomes per nuclei increase with increasing concentration of rapamycin and chloroquine (F) Autophagy spot diameter increases with increasing concentrations of rapamycin and chloroquine.

Starvation is one of the most well-known inducers of autophagy. This study combined chloroquine treatment and serum starvation for two to six hours. Figure 5 demonstrates that increased length of serum starvation results in increased size and number of autophagosomes per nuclei in HeLa cells.

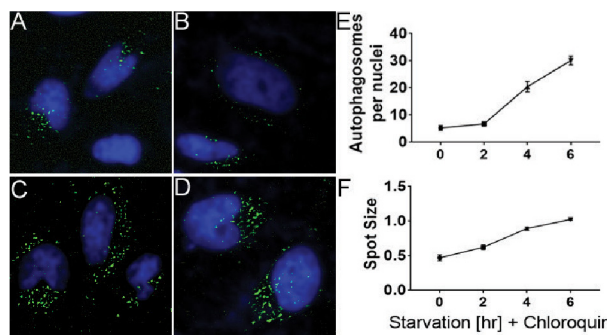


Figure 5. Autophagy-positive spot count increases with longer serum starvation. (A) Control; (B) 2-hour serum starvation + 10 μ M chloroquine; (C) 4-hour serum starvation + 10 μ M chloroquine; (D) 6-hour serum starvation + 10 μ M chloroquine; (E) the number of autophagosomes per nuclei increase according to time in serum-free media. (F) Autophagosome spot diameter increases with increasing time in serum-free media.

Conclusion

The ability to efficiently and rapidly analyze autophagy in living cells is critical for many applications such as screening for compounds that can potentially modify disease states. This application note demonstrates that autophagosome number and size increase in response to known autophagy activator rapamycin and serum starvation. Use of CYTO-ID Autophagy detection kit in combination with Agilent BioTek Gen5 microplate reader and imager software analysis allows consistent and precise measurement of object level data including spot count and spot size.

References

1. Jung, C. *et al.* mTOR Regulation of Autophagy. *Febs Lett.* **2010**, *584*, 1287–1295.
2. Levine, B. *et al.* Autophagy in Immunity and Inflammation. *Nature* **2011**, *469*, 323–35.
3. Shintani, T.; Klionsky, D. J. Autophagy in Health and Disease: a Double-Edged Sword. *Science* **2004**, *306*, 990–5.
4. Coly, P.-M. *et al.* The Autophagy Machinery: A New Player in Chemotactic Cell Migration. *Front. Neurosci.* **2017**, *11*, 78.
5. Henson, E.; Chen, Y.; Gibson, S. EGFR Family Members' Regulation of Autophagy Is at a Crossroads of Cell Survival and Death in Cancer. *Cancers* **2017**, *9*, 27.
6. Scott, R. C.; Schuldiner, O.; Neufeld, T. P. Role and regulation of Starvation-Induced Autophagy in the *Drosophila* Fat Body. *Dev. Cell* **2004**, *7*, 167–78.
7. Levine, B.; Kroemer, G. Autophagy in the Pathogenesis of Disease. *Cell* **2008**, *132*, 27–42.
8. Lorenzi, P. *et al.* A Curated Census of Autophagy-Modulating Proteins and Small Molecules. *Autophagy* **2014**, *10*, 1316–1326.
9. Mizushima, N.; Yoshimori, T.; Levine, B. Methods in Mammalian Autophagy Research. *Cell* **2010**. doi:10.1016/j.cell.2010.01.028
10. Tanida, I. *et al.* Lysosomal Turnover, but not a Cellular Level, of Endogenous LC3 is a Marker for Autophagy. *Autophagy* **2005**. doi:10.4161/auto.1.2.1697
11. Barth, S.; Glick, D.; Macleod, K. F. Autophagy: Assays and Artifacts. *J. Pathol.* **2010**, *221*, 117–24.

www.agilent.com/lifesciences/biotek

For Research Use Only. Not for use in diagnostic procedures.

RA44412.3948611111

This information is subject to change without notice.

© Agilent Technologies, Inc. 2017, 2022
Printed in the USA, March 31, 2022
5994-3378EN