

GFP-LC3 Retroviral Expression Vector

CATALOG NUMBER: RTV-801

STORAGE: -20°C

Components

1. pMX-GFP Control Vector (Part No. RTV-050): One tube, 10 µg at 0.25 µg/µL in TE
2. pMXs-GFP-LC3 Expression Vector (Part No. 380101): One tube, 10 µg at 0.25 µg/µL in TE

Background

Autophagy is a lysosomal degradation pathway for cytoplasmic material, which is activated during stress conditions such as amino acid starvation or viral infection. Mammalian cells use autophagy during short periods of starvation to degrade nonessential cellular components in order to liberate nutrients for vital biosynthetic reactions. Recent results have shown that autophagy also contributes to development, growth regulation and cancer, as well as longevity.

After induction by a stress signal such as amino acid starvation, the first step in autophagy is the formation of an autophagosome. A well published autophagosome marker protein, MAP LC3, was originally identified as a microtubule associated protein and named 'microtubule-associated- protein-light-chain-3'. LC3 is a small 16-18 kDa protein that is soluble in nonstarved cells, but becomes peripherally membrane-associated during amino acid starvation. By immunoelectron microscopy, LC3 has been shown to associate to the inner and outer limiting membranes of autophagosomes, and the membrane association is mediated by a covalent conjugation to a lipid, phosphatidylethanolamine. In Western blots, two forms of LC3 are seen, LC3I and LC3II. LC3I is found in the soluble fraction, and LC3II in the pelletable membrane fraction. Both LC3I and LC3II are seen in nonstarved cells, but during autophagy induction the proportion of LC3II increases. GFP-tagged LC3 expression might be useful as an autophagy assay.

Retroviruses are efficient tools for delivering heritable genes into the genome of dividing cells. Cell Biolabs' pMXs retroviral vector is based on Moloney murine leukemia virus (MMLV). Transfection into a package cell line produces high-titer, replication-incompetent viruses. pMXs-GFP-LC3 is a retroviral expression vector in which human LC-3B gene is fused in frame with GFP and the GFP-LC3 insert is cloned between BamHI and NotI of pMXs vector (Cat.# RTV-010). A pMX-GFP vector without LC3 gene is also provided as a control. The vectors contain the ampicillin-resistance gene, MMLV LTRs, package signal (Figure 1).

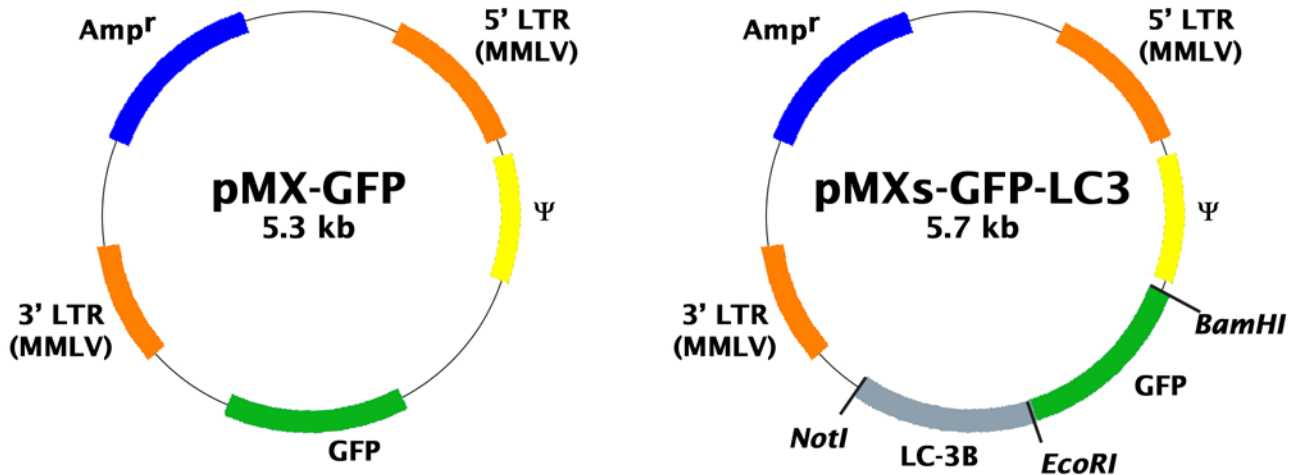


Figure 1. pMX-GFP control vector and pMXs-GFP-LC3 expression vector maps

GFP-LC-3B Sequence in pMXs-GFP-LC3 vector:

```

atggtgagcaagggcgaggagctgttcaccgggggtggtgcccacacctggtcgagctggac
M V S K G E E L F T G V V P I L V E L D
ggcgacgtaaaccggccacaagttcagcgtgtccggcgagggcgagggcgatgccacctac
G D V N G H K F S V S G E G E G D A T Y
ggcaagctgaccctgaagttcatctgcaccaccggcaagctgcccgtgccctggcccacc
G K L T L K F I C T T G K L P V P W P T
ctcgtgaccaccctgacctacggcgtgcagtgttcagccgctaccccgaccacatgaag
L V T T L T Y G V Q C F S R Y P D H M K
cagcacgacttcttcaagtccgccatgcccgaaggctacgtccaggagcgcaccatcttc
Q H D F F K S A M P E G Y V Q E R T I F
ttcaaggacgacggcaactacaagaccgcgagggtgaagttcgagggcgacaccctg
F K D D G N Y K T R A E V K F E G D T L
gtgaaccgcatcgagctgaagggcatcgacttcaaggaggacggcaacatcctggggcac
V N R I E L K G I D F K E D G N I L G H
aagctggagtacaactacaacagccacaacgtctatatcatggccgacaagcagaagaac
K L E Y N Y N S H N V Y I M A D K Q K N
ggcatcaagggtgaacttcaagatccgccacaacatcgaggacggcagcgtgcagctcgcc
G I K V N F K I R H N I E D G S V Q L A
gaccactaccagcagaacacccccatcggcgacggccccgtgctgctgcccgacaaccac
D H Y Q Q N T P I G D G P V L L P D N H
tacctgagcaccagtcggccctgagcaaagacccaacgagaagcgcgatcacatggtc
Y L S T Q S A L S K D P N E K R D H M V
ctgctggagttcgtgaccgccgggatcactctcggcatggacgagctgtacaagtac
L L E F V T A A G I T L G M D E L Y K Y
tcagatctcgagctcaagcttgaattcccatgccgtcggagaagaccttcaagcagcgc
S D L E L K L R I P M P S E K T F K Q R
cgcaccttcaacaaagagtagaagatgtccgacttattcgagagcagcatccaacaaa
R T F E Q R V E D V R L I R E Q H P T K
atcccggtgataatagaacgatacaaggggtgagaagcagcttctgttctggataaaaca
I P V I I E R Y K G E K Q L P V L D K T

```

aagttccttgtagctgaccatgtcaacatgagtgagctcatcaagataattagaaggcgc
K F L V P D H V N M S E L I K I I R R R
ttacagctcaatgctaatacaggccttcttctggtggaacggacacagcatggtcagc
L Q L N A N Q A F F L L V N G H S M V S
gtctccacaccaatctcagaggtgtatgagagtgagaaagatgaagatggattcctgtac
V S T P I S E V Y E S E K D E D G F L Y
atggcttatgcctcccaggagacggttcgggatgaaattgtcagtgtaa
M V Y A S Q E T F G M K L S V -

References

1. Kabeya, Y., Mizushima, N., Ueno, T., Yamamoto, A., Kirisako, T., Noda, T., Kominami, E., Ohsumi, Y., and Yoshimori, T. (2000) *EMBO J.* **19**, 5720-5728.
2. Mizushima, N., Yamamoto, A., Matsui, M., Yoshimori, T., and Ohsumi, Y. (2004) *Mol. Biol. Cell* **15**, 1101-1111.
3. Kabeya, Y., Mizushima, N., Ueno, T., Yamamoto, A., Kirisako, T., Noda, T., Kominami, E., Ohsumi, Y., and Yoshimori, T. (2000) *EMBO J.* **19**, 5720-5728.
4. Mizushima, N., Yamamoto, A., Matsui, M., Yoshimori, T., and Ohsumi, Y. (2004) *Mol. Biol. Cell* **15**, 1101-1111.

License Information

The pMXs vector is licensed from the University of Tokyo.

Warranty

These products are warranted to perform as described in their labeling and in Cell Biolabs literature when used in accordance with their instructions. THERE ARE NO WARRANTIES THAT EXTEND BEYOND THIS EXPRESSED WARRANTY AND CELL BIOLABS DISCLAIMS ANY IMPLIED WARRANTY OF MERCHANTABILITY OR WARRANTY OF FITNESS FOR PARTICULAR PURPOSE. CELL BIOLABS' sole obligation and purchaser's exclusive remedy for breach of this warranty shall be, at the option of CELL BIOLABS, to repair or replace the products. In no event shall CELL BIOLABS be liable for any proximate, incidental or consequential damages in connection with the products.

This product is for RESEARCH USE ONLY; not for use in diagnostic procedures.

Contact Information

Cell Biolabs, Inc.
7758 Arjons Drive
San Diego, CA 92126
Worldwide: +1 858-271-6500
USA Toll-Free: 1-888-CBL-0505
E-mail: tech@cellbiolabs.com
www.cellbiolabs.com

©2009: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.

