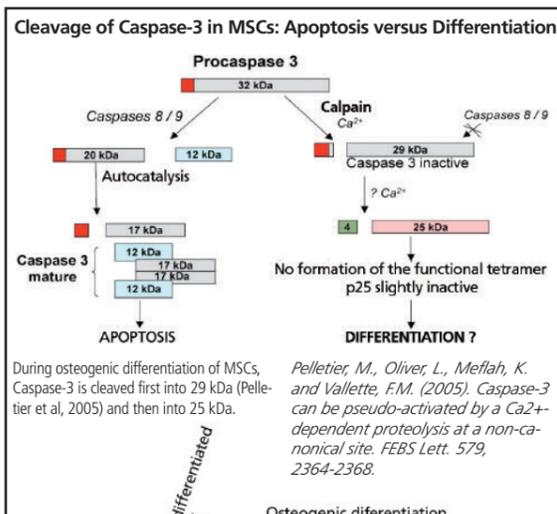


Mesenchymal Stem Cells

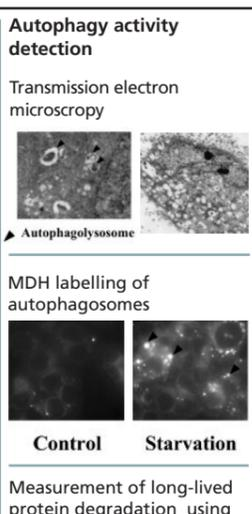
AUTOPHAGY: CELL DEATH OR SURVIVAL? INTERACTION WITH OTHER STRESS PATHWAYS

Oliver L., Hue E., Vallette F. M. Center for Cancer Research Nantes – Angers INSERM UMR 892, University of Nantes, 9 Quai Moncoussu, 44035 Nantes Cedex 01, France

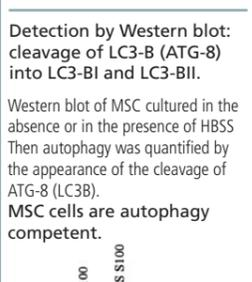
INTRODUCTION
 AUTOPHAGY is a regulated process of the degradation and recycling of cellular constituents, participating in organelle turnover and in the bioenergetics of starvation. Taken to the extreme, autophagy could ultimately result in cell death, through excessive self-digestion and degradation of essential cellular constituents. Thus, it is unclear whether autophagy is fundamentally a cell survival or a cell death pathway – or both.



Autophagy activity detection
 Transmission electron microscopy
 Autophagolysosome
 MDH labelling of autophagosomes
 Control Starvation
 Measurement of long-lived protein degradation using pulse chase ¹⁴C.
 We induce autophagy with HBSS. The % of degraded proteins is compared to the total amount of proteins in cells (addition of 3MA methylamine, an inhibitor for degradation).



Caspase-3 knock-down by shRNA
 Alizarin Red S staining
 Alizarin Red S stains mineral nodules composed of calcium produced by osteoblasts. These are present during terminal osteogenic differentiation. Knock-down of Caspase-3 delays terminal osteogenic differentiation.



The role of Bcl-XL in autophagy in MSCs.
 The essential autophagy protein, Beclin-1 (ATG 6) binds to and is inhibited by Bcl-2 or Bcl-XL. This interaction involves the Bcl-2 homology 3 (BH3) domain in Beclin-1 and the BH3 binding groove of Bcl-2/Bcl-XL.

Effect of Bcl-XL knock-down on autophagic response
 Effect of Bcl-XL knock-down (note that little or no Bcl-2 is detectable in MSCs) MSCs were infected with shBcl-XL or src viral particles. 48 h later total cell extracts were made and Western Blot determined the expression of Bcl-XL. The efficiency of the knockdown of Bcl-XL was about 80% (average of 3 experiments) as compared to an internal control actin.

Alizarin Red S staining
 RT-qPCR analysis of osteogenic markers during differentiation.
 MSCs are not differentiated, there are no markers for osteoblasts. If we block Caspase-3 this differentiation is disrupted. If we add growth factors and block Caspase-3, the cells try other ways to induce differentiation: the genes for osteogenic differentiation RUNX2, OPN and MGP are activated. (indirect results)

MDH staining of MSCs
 MSC-Bcl-XL cells and MSC-scr cells were cultured in HBSS for 6 h to induce autophagy in these cells. The cells were cultured for the last 30 min in the presence of 1 μM Monodansyl pentane (MDH) to label autophagosomes. At the end of the incubation, the cells were washed with PBS then analysed under a microscope at Ex: 335 nm/Em: 525 nm.

Effect of Bcl-XL knock-down on autophagic response Confirmation of the results.
 Since the knock-down in MSC cells affects between 60-80% of the cells, we need to confirm the results. Therefore knock-down was done using a pSilencer shRNA GFP-Bcl-XL. After transfection the cells were cultured in the presence or in the absence of HBSS for 6 h and the stained with MDH.

Graph of percentage knock-down of Bcl-XL.
 MSCs were cultured in the presence of HBSS and the cells were photographed every 10 min over 72 h. The number of dead cells was quantified at each point. The graphs depict the results from 3 different experiments: 3 different cells of MSCs. We see here that autophagy is not accompanied by cell death.

Conclusion
 In MSCs, surprisingly Bcl-XL stimulated the autophagic activity in cells, which could be qualified by an increase in the number and size of the autophagosomes. The knock-down of Bcl-XL completely inhibited autophagic activity and sensitized these cells to apoptotic insults. Finally, in MSCs, autophagy appears to be the predominant stress sensor mechanism, and is likely to be used as a protection against apoptotic insults.