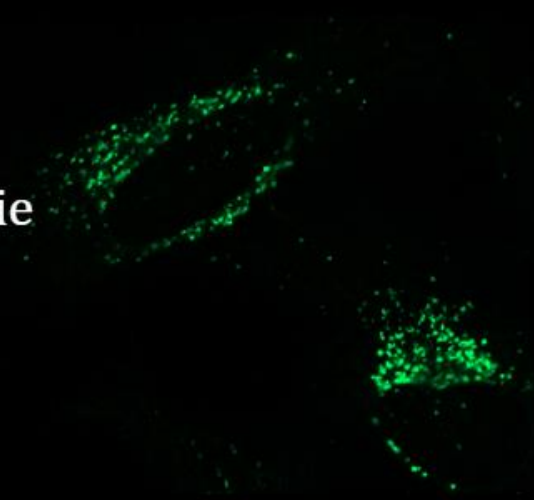
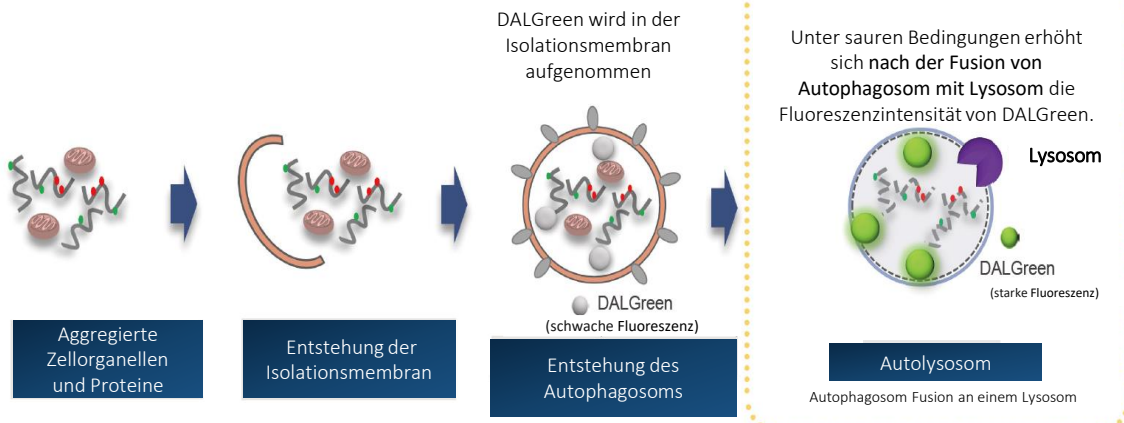


# Visualisierung von Autophagie

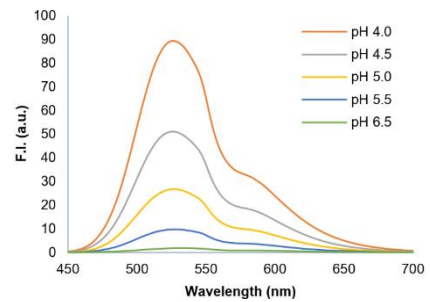
D675 DALGreen - Autophagy Detection



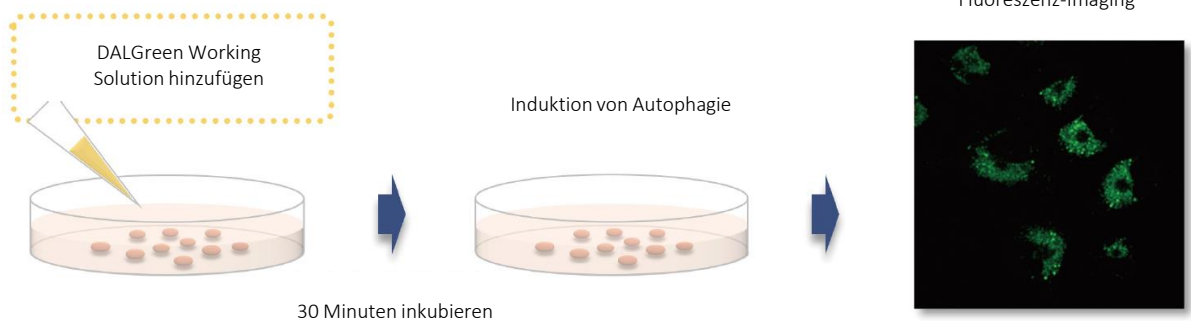
## ✓ Neue Technologie



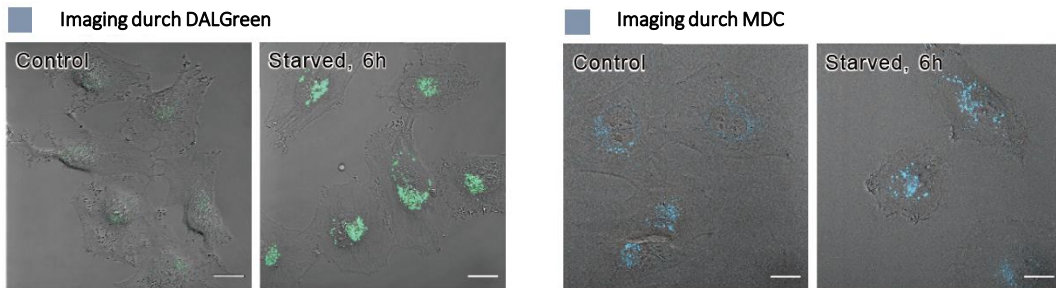
DALGreen ist ein pH-abhängiger Farbstoff (starke Fluoreszenzintensität ca. pH 4,5)



## ✓ Einfaches Verfahren



## ✓ Besseres Signal

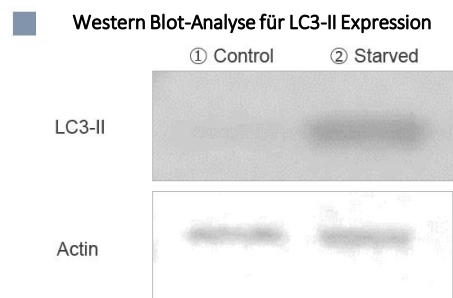
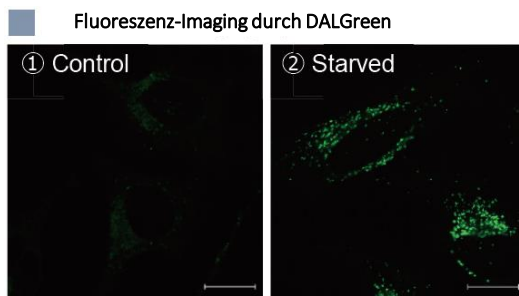


### Quelltext : 3<sup>rd</sup>. Autophagy Guideline

pharmacologically up- or downregulated.<sup>991,992</sup> Whether this method can also be used in adult animals needs to be determined. Furthermore, it should be kept in mind that staining with MDC is not, by itself, a sufficient method for monitoring autophagy (see Acidotropic dyes).

As stated above with regard to other organisms, staining with MDC or derivatives (such as monodansylmethylamine) is not sufficient for detection of autophagy, as these stains also detect vacuoles. The same is true for the use of LysoTracker Red, Neutral Red or acridine orange. The fluorophore of the

## ✓ Gute Korrelation mit LC3



### DAL Green – Autophagy Detection

Code #            Packungsgröße  
D675-10        20 nmol



Für weitere Informationen besuchen Sie unsere Website <http://www.dojindo.eu.com/store/p/913-DALGreen-Autophagy-Detection.aspx>

**DOJINDO** STATE OF THE ART  
LIFE SCIENCE TECHNOLOGIES

EUROPEAN HEADQUARTERS  
DOJINDO EU GMBH  
Leopoldstr. 254, 80807 Munich, Germany  
Phone            +49 89 3540-4805  
Fax                +49 89 3540-4806  
email             [info@dojindo.eu.com](mailto:info@dojindo.eu.com)  
[www.dojindo.eu.com](http://www.dojindo.eu.com)

DISTRIBUTED BY