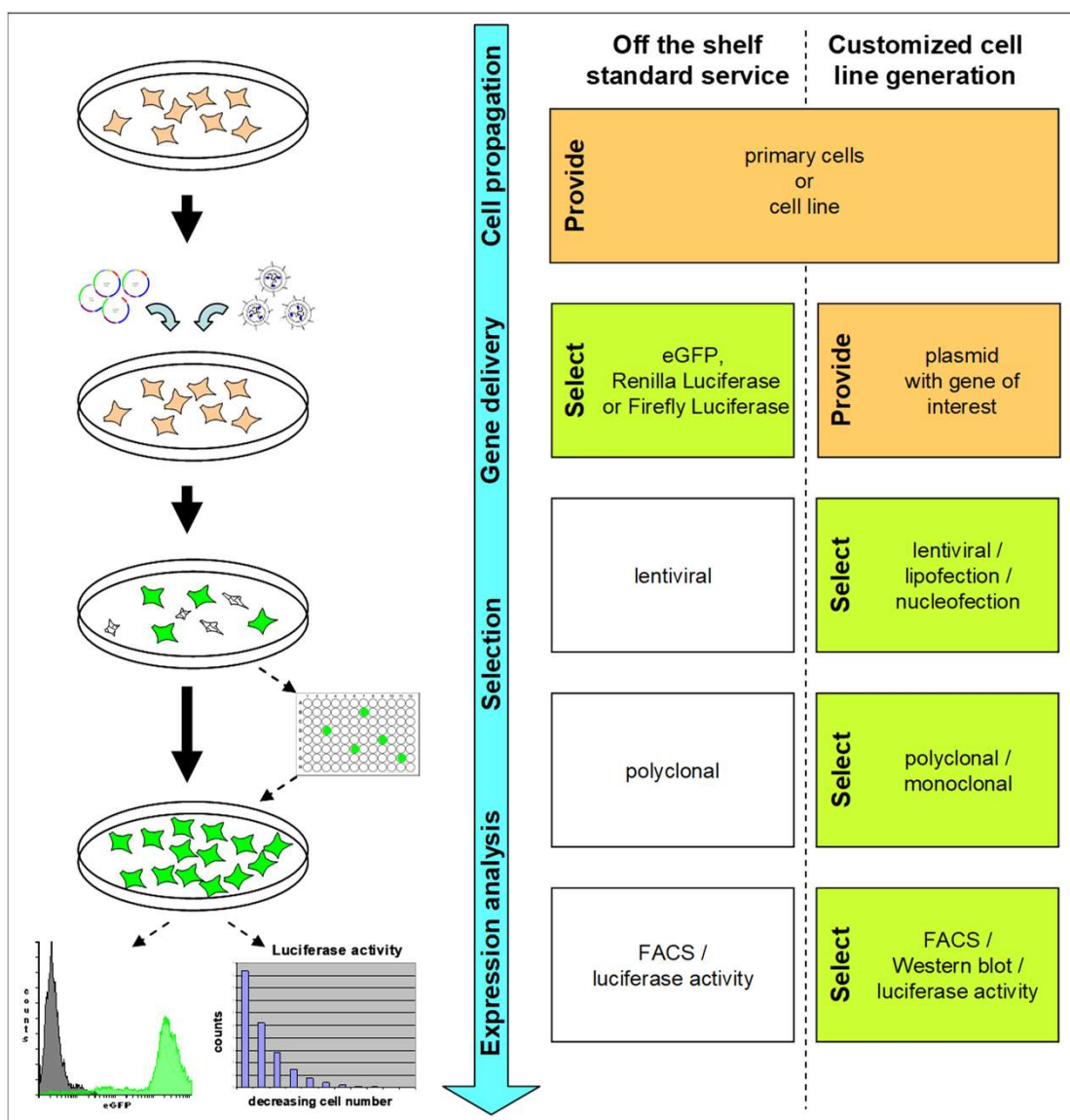


## ➤ Stable generation of cell lines and primary cells

Stable generation of cell lines can serve many purposes, either looking for the functional impact of the knock-in or knock-out of genes of interest, or just for labeling the cells. Especially the labeling of cell lines becomes more important in order to track cells in vivo or to perform proliferation studies based on luciferase-expression. Based on extensive hands-on experience in genetically engineering cells, the „Stable cell line generation service“ provides the possibility for genetic modifications using a.) standard transfection methods (lipofection/nucleofection/Ca2-PO4-transfection) or b.) lenti-viral transduction which allows for highly efficient genetic manipulation not only of cell lines but also of primary cells and non-dividing cells.

## ➤ Cell Line Generation Procedure



Contact us to discuss your project: [info@proqinase.com](mailto:info@proqinase.com)

ProQinase disclaims any warranty explicitly or implied that the use of this service is free from third party intellectual property claims unless this is explicitly stated.