



# ID-Centrifuge L ■ ID-Incubator L

The Next Generation

