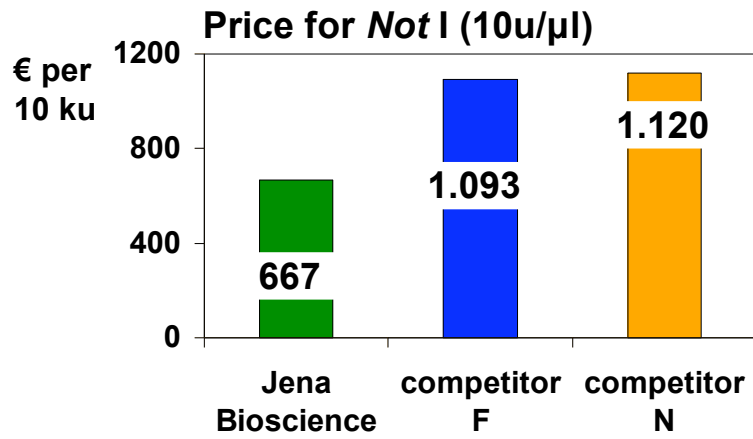
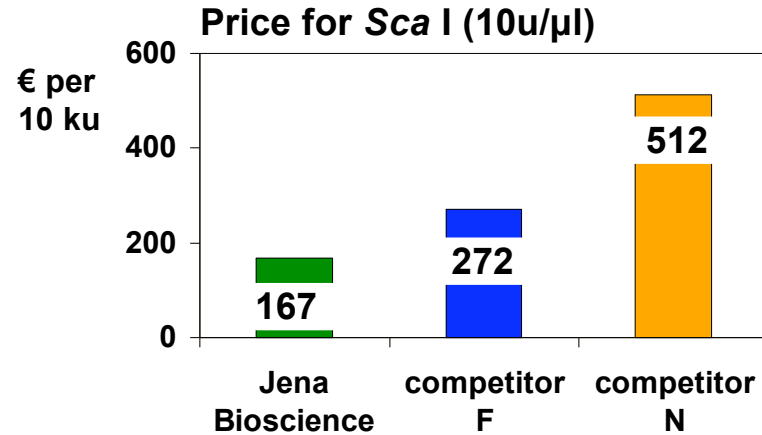


10.000 units of restriction enzyme can cost a fortune or be a bargain*:



Save > 40% with
Jena Bioscience enzymes



Save > 67% with
Jena Bioscience enzymes

Attractive discounts on our list prices available, just contact
enzymes@jenabioscience.com for details

* prices based on: Jena Bioscience list prices, Competitor F prices according to pricelist for Germany valid since May 15, 2006; Competitor N prices taken from competitor N's German web site