

Protocol

Quick & Easy Yeast Transformation Mix



The Quick & Easy Yeast Transformation Mix (Cat. No. 631851) Visual Protocol is provided for your convenience, but is not intended for first-time users. For additional details, see the Quick & Easy Yeast Transformation Mix Protocol-At-A-Glance.



1

Pick a colony or colony-sized amount of yeast from a culture plate



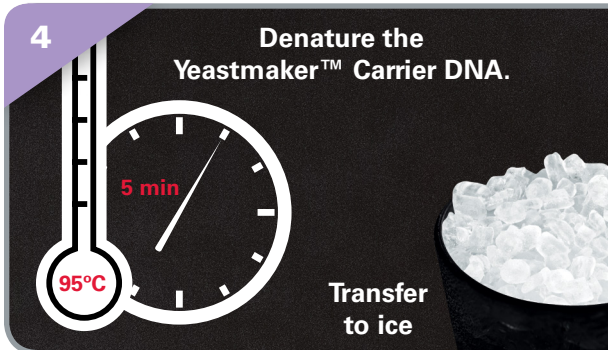
2

Resuspend yeast cells in 300 µl sterile water



3

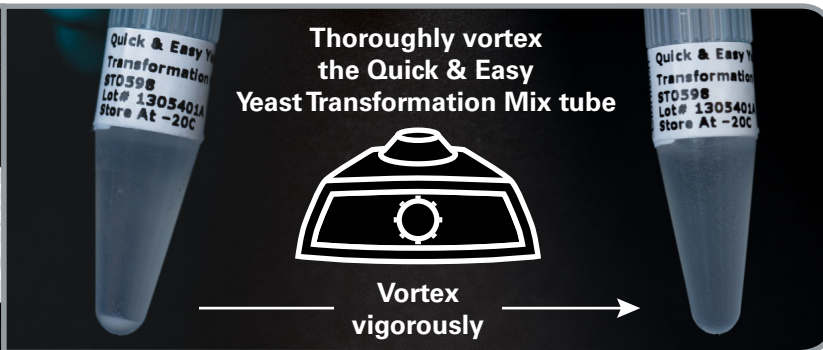
Centrifuge the cells at 3,000 x g for 3 min, and carefully remove supernatant



4

Denature the Yeastmaker™ Carrier DNA.

Transfer to ice



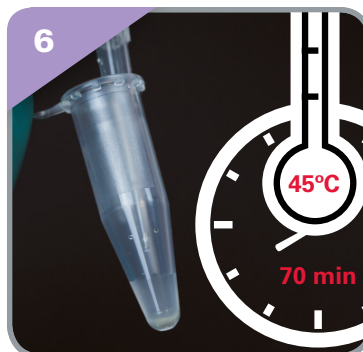
Thoroughly vortex the Quick & Easy Yeast Transformation Mix tube

Vortex vigorously

Prepare individual reagents (can be done in advance)

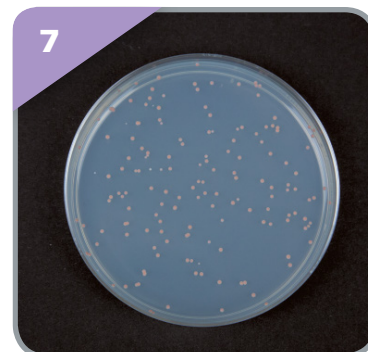
5 **Prepare transformation mixture**

Vol/Rxn	# Rxns	=	Reagent
93.5–94 µl	x _____	= _____	Quick & Easy Yeast Transformation Mix
5 µl	x _____	= _____	Yeastmaker Carrier DNA (Denatured)
1.0–1.5 µl	x _____	= _____	Transforming DNA (150–200 ng)
100 µl	x _____	= _____	Total Volume



6

Resuspend pellet in 100 µl transformation mixture, and incubate



7

Dilute the transformations 10-fold and 100-fold with sterile water, and plate