

RapidTrans™

efficient, affordable competent *E. coli* in a convenient format

RapidTrans™ are high-efficiency competent *E. coli* supplied in a convenient 96-tube tray. Ideal for cloning, plasmid preparation and library construction, RapidTrans cells are affordably priced and packaged in a format that provides maximum flexibility while eliminating waste. Each tube contains 50 µl of cells for one transformation reaction. In contrast to other 96-well formats that require thawing all 96 reactions at one time, RapidTrans enables the use of as few or as many reactions as needed, without thawing the other cells. This eliminates the reduced efficiencies and waste caused by repeated freeze/thaw cycles.

The RapidTrans™ Kit advantage

- Consistent high-quality, high-efficiency *E. coli*
- Freedom to use anywhere from 1-96 reactions eliminates wasted reactions
- Economical pricing
- Adaptable to high-throughput use

Strains and Genotypes

TAM1: *mcrA* Δ (*mrr-hsdRMS-mcrBC*) Φ 80*lacZ* Δ M15
 Δ *lacX74 recA1 araD139 Δ (*ara-leu*)7697 *galU galK*
*rpsL endA1 nupG**

TAM1 *E. coli* have been genetically modified for use in cloning, plasmid preparation and library construction. They are offered either as Competent *E. coli* at $> 1 \times 10^8$ cfu/µg.



Product	Format	Catalog No.
RapidTrans™ TAM1	1 x 96 rxns	11096
Competent <i>E. coli</i>	5 x 96 rxns	196

CONTENTS & STORAGE

One or five 96-tube trays (50 µl/tube) of competent *E. coli*, pUC19 plasmid DNA and SOC medium. Storage conditions vary from room temperature to -80°C, see manual for details.