

# Transformation Experts

Chemi-Competent Cells



Introducing the emerging standard in competent cells  
from Bio-Rad, the transformation experts.

**BIO-RAD**

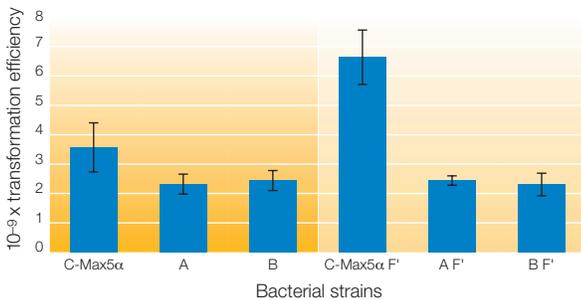


# Chemi-competent cells for high-efficiency transformation

## C-Max™5α and C-Max5α F' Chemi-Competent Cells

C-Max5α and C-Max5α F' *E. coli* strains are provided in convenient 0.2 ml or 0.05 ml single-use aliquots to maintain consistent performance without loss of transformation efficiency due to repeated freeze-thawing. Simply thaw the cells, add DNA, and complete your protocol.

Both cell types are perfect for many cloning applications, including those requiring large plasmid constructs. C-Max5α F' chemi-competent cells provide all of the features of C-Max5α cells but also contain an F' episome and the *lacI<sup>q</sup>* gene, which permits regulation of *lac*, *tac*, and *trc* promoters.



Comparison of transformation efficiency of chemi-competent cells. A and B are comparable strains from the top two suppliers of competent cells.

## C-Max5α Chemi-Competent Cells

**Genotype:** F<sup>-</sup> φ80dlacZΔM15 Δ(lacZYA-argF)U169 recA1 endA1 hsdR17 (r<sub>k</sub><sup>-</sup>, m<sub>k</sub><sup>+</sup>) phoA supE44 λ<sup>-</sup> thi-1 gyrA96 relA1

- Transformation efficiencies >1 × 10<sup>9</sup>
- Mutations facilitating transfection with large plasmids
- Blue/white screening for recombinants
- *endA1* and *recA* mutations for high quality, high plasmid copy numbers, and a lowered risk of recombination

## C-Max5α F' Chemi-Competent Cells

**Genotype:** φ80dlacZΔM15 Δ(lacZYA-argF)U169 recA1 endA1 hsdR17 (r<sub>k</sub><sup>-</sup>, m<sub>k</sub><sup>+</sup>) phoA supE44 λ<sup>-</sup> thi-1 gyrA96 relA1/F'[[*lacI<sup>q</sup>* Tn10 (Tet<sup>r</sup>)]

- Transformation efficiencies >1 × 10<sup>9</sup>
- Mutations facilitating transfection with large plasmids
- Regulatable expression due to the *lacI<sup>q</sup>* gene
- Blue/white screening for recombinants
- *endA1* and *recA* mutations for high quality, high plasmid copy numbers, and a lowered risk of recombination
- Susceptible to infection by bacteriophage M13

## Ordering Information

Catalog #	Description
170-3340	C-Max5α Chemi-Competent Cells, 5 x 0.2 ml
170-3342	C-Max5α Chemi-Competent Cells, 10 x 0.05 ml
170-3341	C-Max5α F' Chemi-Competent Cells, 5 x 0.2 ml
170-3343	C-Max5α F' Chemi-Competent Cells, 10 x 0.05 ml



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** www.bio-rad.com **USA** (800) 4BIORAD **Australia** 02 9914 2800 **Austria** (01)-877 89 01 **Belgium** 09-385 55 11 **Brazil** 55 21 507 6191  
**Canada** (905) 712-2771 **Czech Republic** + 420 2 41 43 05 32 **China** (86-21) 63052255 **Denmark** 44 52 10 00 **Finland** 09 804 22 00  
**France** 01 47 95 69 65 **Germany** 089 318 84-177 **Hong Kong** 852-2789-3300 **India** (91-124)-6398112/113/114, 6450092/93 **Israel** 03 951 4127  
**Italy** 39 02 216091 **Japan** 03-5811-6270 **Korea** 82-2-3473-4460 **Latin America** 305-894-5950 **Mexico** 52 5 534 2552 to 54  
**The Netherlands** 0318-540666 **New Zealand** 64 9 415 2280 **Norway** 23 38 41 30 **Poland** + 48 22 8126 672 **Portugal** 351-21-472-7700  
**Russia** 7 095 721 1404 **Singapore** 65-62729877 **South Africa** 00 27 11 4428508 **Spain** 34 91 590 5200 **Sweden** 08 555 12700  
**Switzerland** 061 717-9555 **Taiwan** (8862) 2578-7189/2578-7241 **United Kingdom** 020 8328 2000