



Transformation Experts

Introducing the emerging standard in
high-efficiency competent cells
from Bio-Rad, the leader
in gene transfer technology.



BIO-RAD

Competent cells for expert transformation



- High efficiencies
- Consistent performance
- Blue/white screening for recombinants
- Optimized volumes

Chemi-Competent Cells

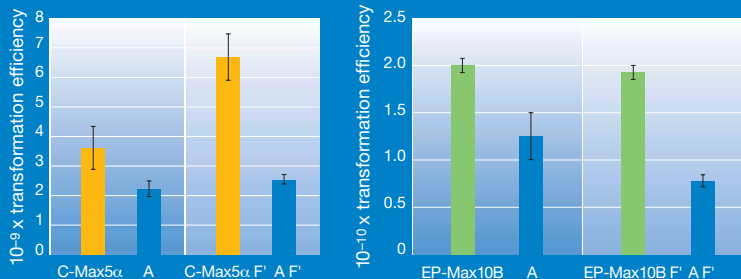
C-Max™ 5α (>1 x 10⁹/μg DNA)

C-Max5α F' (>1 x 10⁹/μg DNA)

Electro-Competent Cells

EP-Max™ 10B (>1 x 10¹⁰/μg DNA)

EP-Max10B F' (>1 x 10¹⁰/μg DNA)



Comparison of transformation efficiency with comparable strains from a leading supplier of competent cells. Cells were transformed with pUC19 DNA.

BIO-RAD