

## Statement of nonpathogenicity

Product	Cat. No.
ccdB Survival 2 T1 <sup>R</sup> cells	A10460
DH5α cells	12034-013, 18288-019, 18263-012, 11319-019, 12297-016, 18258-012, 18265-017
DH10B cells	C5100-03, 18290-015, 18297-010, 12033-015, 12331-013
OmniMAX 2 T1 <sup>R</sup> cells	C8540-03
TOP10 cells	C4040-10, C4040-03, C4040-06, C4040-52, C664-11, C664-55, C664-24
TOP10F' cells	C3030-03, C3030-06, C665-11, C665-24
BL21 cells	C6010-03, C6020-03, C6070-03, C6565-03, C6060-10, C6060-03, C6000-03

Invitrogen™ ccdB Survival™ 2 T1<sup>R</sup>, DH5α™, DH10B™, OmniMAX™ 2 T1<sup>R</sup>, TOP10, and TOP10F' competent cells are derived from *E. coli* K12. K12-derived strains are unable to colonize and are therefore considered to be nonpathogenic in humans or other animals. For further information on this topic, please review the K12 safety literature such as that found at [epa.gov/sites/production/files/2015-09/documents/fra004.pdf](http://epa.gov/sites/production/files/2015-09/documents/fra004.pdf)

BL21 cells are derived from a B strain of *E. coli* and are considered to be nonpathogenic. While the potential for colonization is slightly greater than that of K12 strains, they remain a Risk Group 1 agent and should be handled using Biosafety Level 1 conditions.

**Carla Potts**

Product Manager, Competent Cells  
Thermo Fisher Scientific

Find out more at [thermofisher.com/compcells](http://thermofisher.com/compcells)

