

## DIMENSIONS OF CCS LID WITHOUT SPACERS

*PN 6000025 --- clear, non-sterile lid*

*PN 6000024 --- clear, sterile lid*

*PN 6000027 --- black, non-sterile lid*

*PN 6000026 --- black, sterile lid*

### Material and treatment

Black or clear polystyrene CCS lids without stacking spacers. Lids are available sterile or non-sterile. Universal design (no corner cut-offs). Lids are developed for PerkinElmer's miniTrak, plateTrak, or PlateStak applications.

### Packaging

Lids are packed in cases of 200 pieces, packed as 8 bags of 25 lids each.

### Dimensions



Outside view black CCS lid without spacers



Inside view black CCS lid without spacers

Lid Dimensions	6000024/-25/-26/-27
Inside length	125.3 mm
Inside width	83.1 mm
Maximum height	9 mm