



## 18 mm Circular Cover slides – plain

circle diameter: 18 mm

thickness: 0,22 mm

Cat. No.	Amount
CSL-103	1.000 pcs.

**For general laboratory use.**

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at ambient temperature

**Shelf Life:** n/a