



## 22 mm Thick Circular Cover slides - siliconized

circle diameter: 22 mm

thickness: 0,5 mm

Cat. No.	Amount
CSL-107	500 pcs.

**For general laboratory use.**

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at ambient temperature

**Shelf Life:** n/a