



22 mm Circular Cover slides – siliconized

circle diameter: 22 mm

thickness: 0,22 mm

Cat. No.	Amount
CSL-106	500 pcs.

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: n/a