



## Puck Support

SPINE

Cat. No.	Amount
CC-CSM003-0005A	1 pc.

**For general laboratory use.**

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at ambient temperature

**Shelf Life:** n/a