

PRODUCT NAME 2 ml Microtubes for SX-8G IP-Star® Compact		
Cat. No: WA-008-0100	Lot #: 001	Format: 100 tubes
Remark: Low Binding		

Description

These tubes are intended for being used on the SX-8G IP-Star® Compact Automated System for dispensing the reagents during the corresponding automated protocols.

Characteristics

- Volume: 2 ml
- Diameter: 10.8 mm
- Length: 46 mm
- Base shape conical
- Type of closure screw cap
- Material: polypropylene [PP]
- Colour: transparent
- Non-sterile tube
- Screw cap enclosed