



sartorius stedim
biotech

Accessories for Adapter Sanitary 1ZA0GV0003

EPDM rings for 3/4" Tri-Clamp, 2 items
Order No.: 17521 - - - 013

Tri-Clamp, 1 item
Order No.: 7ZSB - - - 0001