

# Accessories for Adapter Sanitary 1ZAOGV0003

EPDM rings for  $\frac{3}{4}$ " Tri-Clamp, 2 items  
Order No.: 17521---013

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