

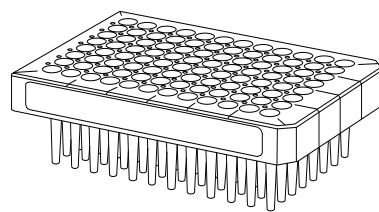
SOLA μ SPE well plates

Positive pressure manifold set-up considerations

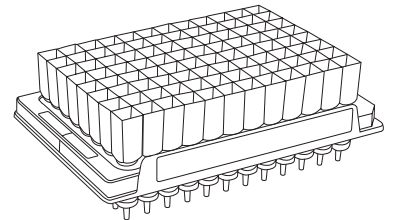


1

The Thermo Scientific™ SOLA μ ™ solid phase extraction (SPE) well plate is used in exactly the same manner as standard micro elution plates. However, not all micro elution SPE plate formats are the same. Some adjustment of the manifold maybe required.



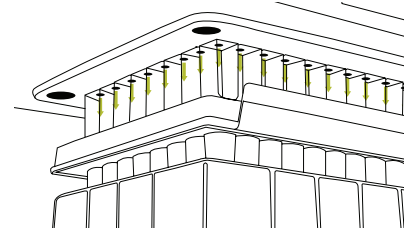
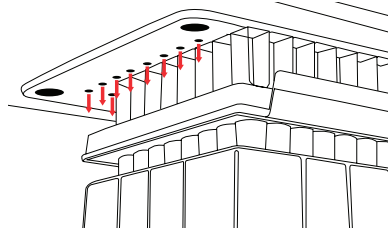
Competitor μ well plate



SOLA μ SPE well plate

2

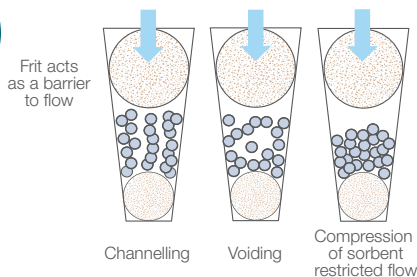
- Make sure the SPE plate fits directly onto the collection plate
- Make sure the tips of the SPE plate fit into the wells of the collection plate
- Make sure the Positive Pressure seal is correctly aligned with the top of the SPE plate
- Prior to starting the SPE process make sure the pressure is off
- Seal the plate and increase the pressure until flow is observed



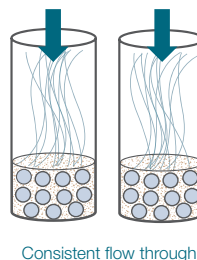
- **Note** – you may experience more controlled flow and a requirement for higher pressure with the SOLA μ plate (due to its advanced fritless design) which allows for better interaction with the sorbent and sample

3

Competitor μ well plate



SOLA μ SPE well plate



Flow-through characteristics: SOLA μ vs. Competitor μ well plates

- Consistent manufacture ensures the same amount of material is packed every time
- Frits act as a barrier to flow, removal aids flow
- Space between media is not controlled by packing density
- SOLA unit structure has increased porosity due to manufacture control

Find out more at thermofisher.com/solaspe