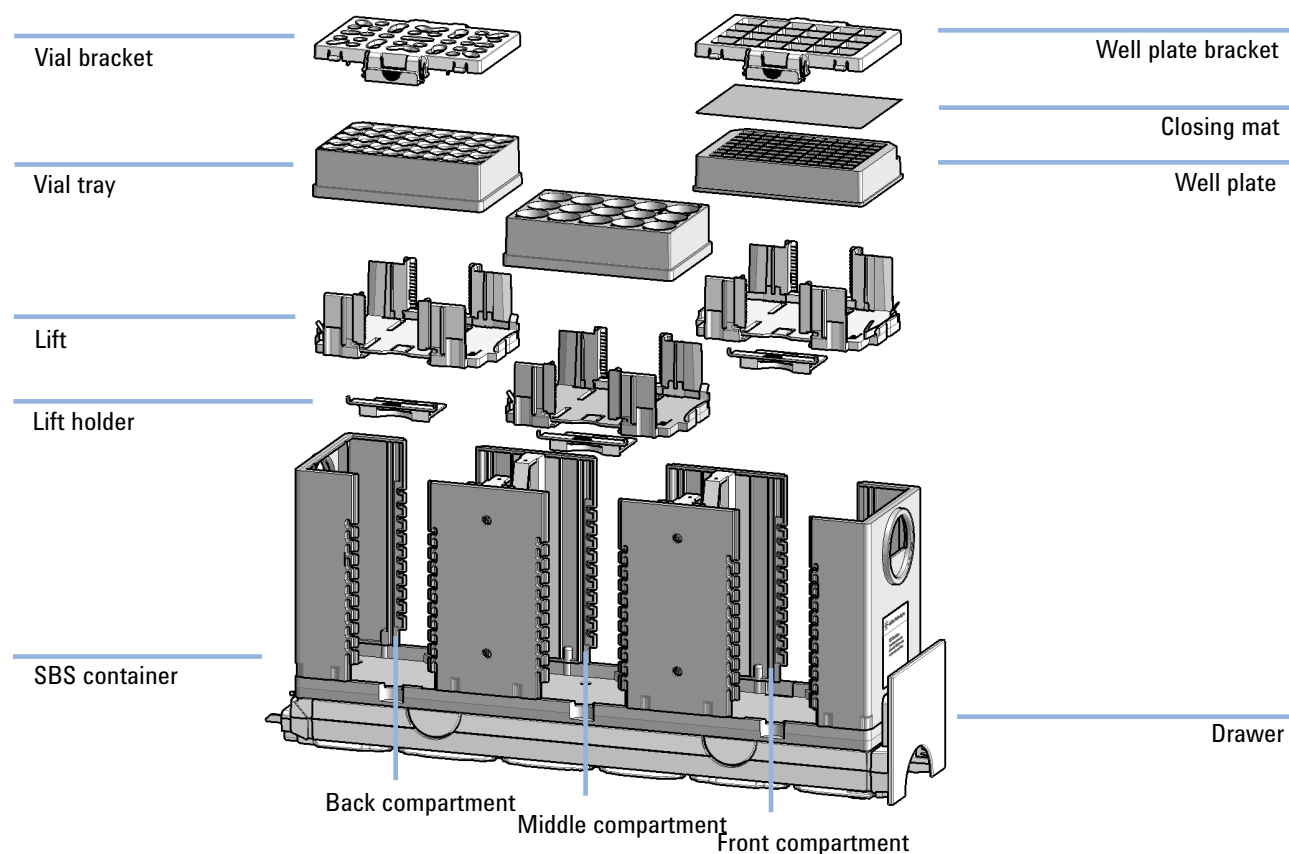


# Agilent 1290 Infinity II Preparative Open-Bed Sampler/Collector SBS Container Configuration and Application

## Technical Note

This technical note describes the configuration and application options of the SBS container for the Agilent 1290 Infinity II Preparative Open-Bed Sampler/Collector (G7158B) and Agilent 1290 Infinity II Preparative Open-Bed Fraction Collector (G7159B). This includes operation and modification of the container, usage of brackets for sampling and correct height determination of the lifts.

## Configure the SBS Container



**Figure 1** Overview of SBS container and accessories



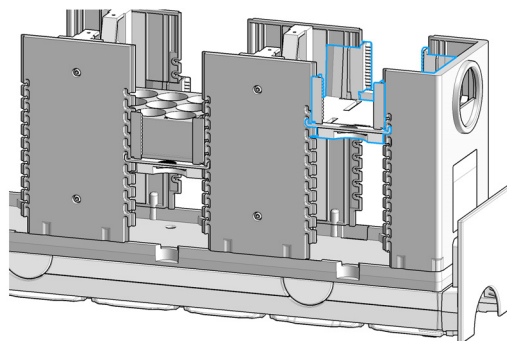
- 1 After transport, the lift and the lift holders are at the bottom of the container compartment. Pull the lift up until the lift holders are accessible. Move the lift holders up until they lock in the lowest position.

### CAUTION

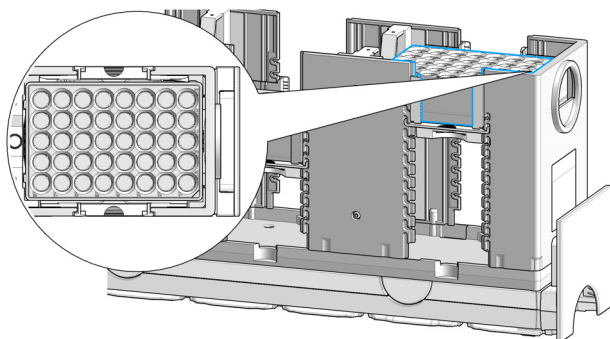
#### Protruding objects

Objects (for example, vials, tubes, brackets) protruding over the upper edge of the SBS container can damage or brake the injection needle.

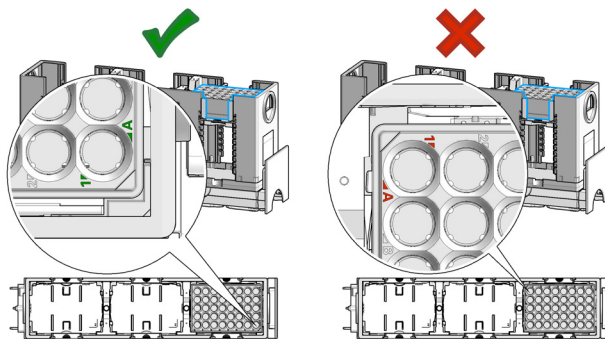
- Mount all sample containers at a level where they do not protrude beyond the upper edge of the SBS container. Take seals, caps, and brackets into consideration.
  - Agilent Technologies recommends that you use Agilent products for which the level in the compartment is specified.
- 2 Pull out the lift holders, move the lift upward, and push the lift holders in again to align the lift at the desired height. For a detailed description of the sample container types and their position levels in the compartment, see [Table 1](#) on page 4.



- 3 Place the sample container (for example, vial tray, well plate or sample holder) in the lift.



- 4 Ensure that the sample containers in the compartments are correctly positioned and oriented (A1 on A1).



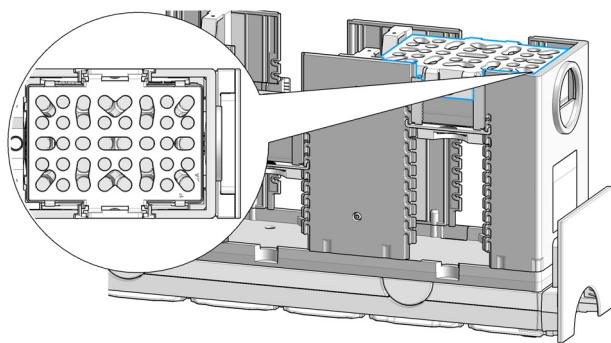
### CAUTION

#### Pinched needle

The injection needle can get trapped in the cap of sample vials or in the mat of well plates.

→ Cover the vial trays and well plates with the appropriate bracket.

5 Prepare the vial tray and cover it with the corresponding bracket.



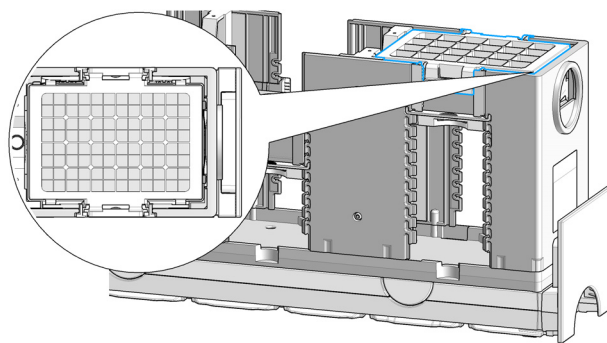
### CAUTION

#### Pinched needle

The injection needle can get trapped in the cap of sample vials or in the mat of well plates.

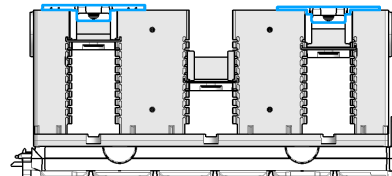
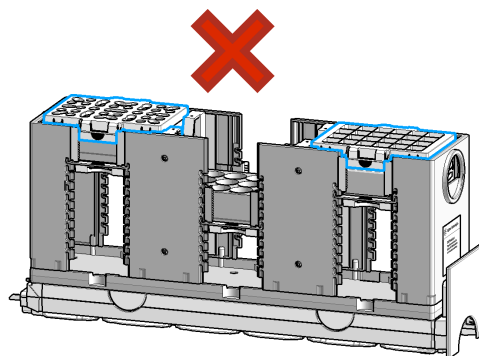
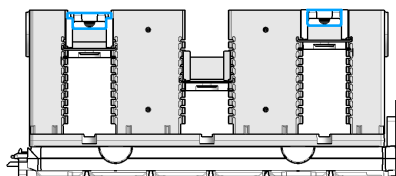
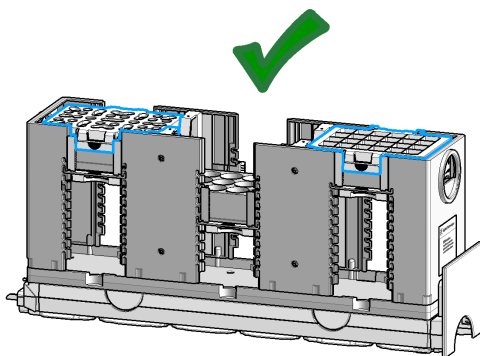
→ Cover the vial trays and well plates with the appropriate bracket.

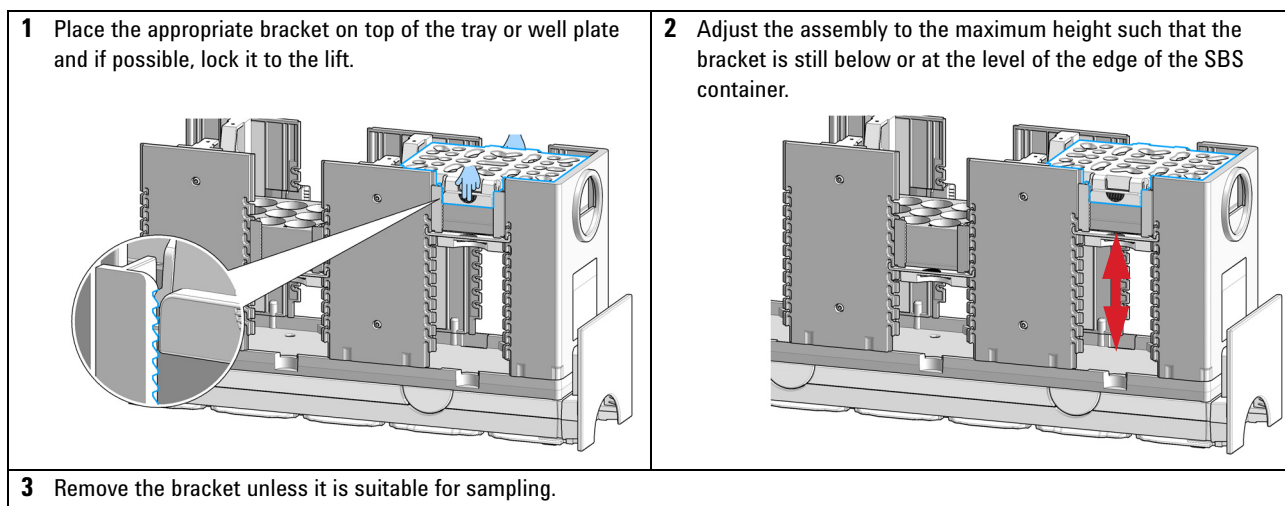
6 Cover the well plate with the mat and the corresponding bracket.



## Determine the Correct Level of the SBS Lift

For determination of the correct level of the SBS lift, always employ either the well plate bracket for well plates or the vial bracket for all other kinds of trays. Sample containers must not protrude beyond the upper edge of the SBS container.





The required levels for Agilent supported trays and well plates are listed in [Table 1](#) on page 4.

## SBS Container Plates

**Table 1** Overview of SBS container plates

Agilent part number	ChemStation name	Name	Compatible sealing	Max. allowed fill volume [mL]	SBS level
5043-9300	A_5043-9360_96_2.2mL	96 wellplate 2.2 mL		1.98	2nd highest
5043-9305	A_5043-9362_96_1.0mL	96 wellplate 1.0 mL		0.90	highest
5043-9308 5043-9309	A_5043-9363_96_1.2mL	96 wellplate 1.2 mL		1.00	highest
5023-2471	40VialPlate	2 mL Vial Plate	5185-5824	1.50	2nd highest
5043-1826	5043-1826 15x6mL Vials	5 mL HR and 6 mL Vial Plate	5188-2758 9301-1379	5.00 6.00	2nd highest

### NOTE

For features such as dilution and mixing, adjust fill volumes accordingly if the sample tends to foam.

