

Mounting Guideline

Sirius L standard sample holder (for 75x12mm and 55x12 mm tubes)

The standard sample holder (for 75x12mm and 55x12mm tubes) is preinstalled in the factory for use with 75mm tubes. For use with 55mm tubes follow the instructions below. Please refer to the user's manual chapter 2.1 and 3.4.1 for further information, too.

A: For 75mm tubes:

- use the preinstalled standard sample holder only.

B: When using 55 mm tubes:

- unscrew the standard sample holder
- add the clear cylinder from the bottom to the reflector
- place back and screw the reflector with insert into the instrument!

