

trUView Beam Height Adaptor

To install adaptor:

1. Screw beam height adaptor on the installation tool provided.
2. Lower beam height adaptor into cuvette holder.
3. Unscrew installation tool.

To remove adaptor:

1. Screw installation tool onto beam height adaptor.
2. Lift adaptor out.

