

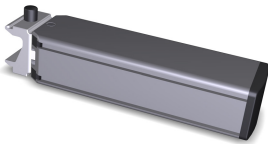
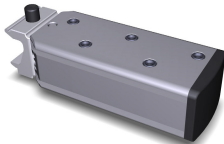
# Attach rail extension and column holders




## Instructions for Use

### 1 Scope

Instructions on how to attach the following rail extension and column holders to the instrument.

**Note:** *The described column holders are designed for the ÄKTA™ avant, ÄKTA pure™, and ÄKTA go™ instruments and may not be possible to attach to other ÄKTA instruments. Also, column holders used for other ÄKTA instruments may not be possible to attach to the ÄKTA avant, ÄKTA pure, and ÄKTA go instruments.*

Holder	Illustration
Rail extension (29011352)	
Column holder rod (28956270)	

Holder	Illustration
Column holder (28956282)	
Column clamp o.d. 10-21 mm (28956319)	
Multi-directional column clamp (29383530)	

## 2 Attach Rail extension and Column holder rod

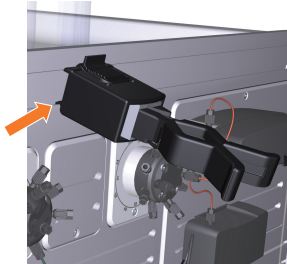
Press down the button on the rod (1) and attach the rod to the rails of the instrument (2).



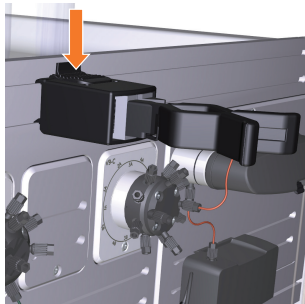
### 3 Attach Column holder

Step	Action
------	--------

- 1 Attach the lower hooks of the column holder snap-in to the lower rail.



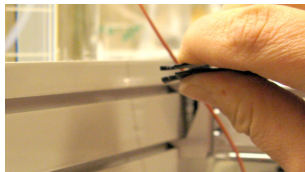
- 2 Press the tab on the holder and attach the upper hooks of the column holder snap-in to the upper rail.



### 4 Attach Column clamp

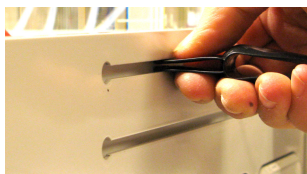
Step	Action
------	--------

- 1 Press the inner end tabs of the clamp together.

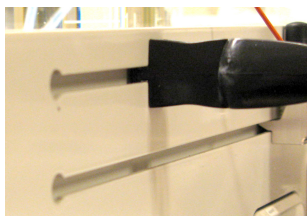


Step	Action
------	--------

- |   |   |
|---|---|
| 2 | Insert the clamp into a rail of the instrument. |
|---|---|



- |   |  |
|---|--|
| 3 | Turn the clamp 90 degrees and release. |
|---|--|

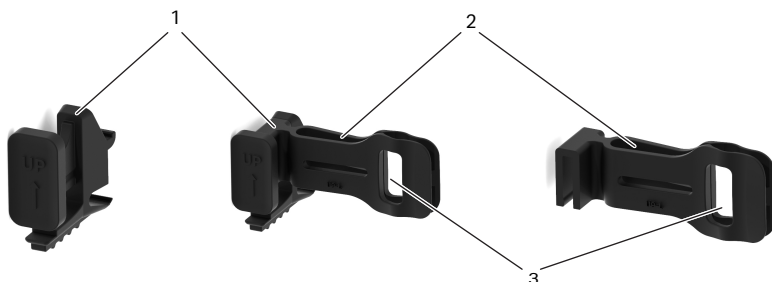


## 5 Multi-directional column clamp

The Multi-directional column clamp can be used to attach small sized columns with an outer diameter of 12 to 18 mm directly to the UV monitor. The direction of the clamp can be changed for holding columns directly to UV monitor **U9-M**, **U9-T**, or **U9-L**.

Two base units of the Multi-directional column clamp can be used to mount a column horizontally on instruments with long rails.

The illustration below shows the Multi-directional column clamp.



Part	Description
1	Base unit with snap-in to holder rails
2	Clamp unit
3	Position for column

**Tip:** Orient the clamp so that the name of the UV monitor, **U9-M** or **U9-L**, can be read on the front of the clamp.

- For **U9-M**, place the base clamp to the right of the column. Make sure that the text **U9-M** shows on the front of the clamp.
- For **U9-T**, and **U9-L**, place the base clamp to the left of the column. Make sure that the text **U9-L** shows on the front of the clamp.

## 5.1 Mount the column

Step	Action
------	--------

- |   |  |
|---|--|
| 1 | Attach the base unit: <ol style="list-style-type: none"><li>Press the upper tab into the upper rail.</li></ol> |
|---|--|

Step	Action
------	--------

- b.** Press the lower tab towards the lower rail to insert it.



*Result:*



**Note:**

*To remove the base unit, press the bottom tab upwards and take the base unit out from the rails.*

- 2 To mount the column horizontally:
  - a.** Attach a second base unit to the instrument.
  - b.** Place the column on the two base units.

Step	Action
------	--------

c. Adjust the location of the base units to support the column.





3 To mount the column vertically, attach the clamp unit on the base unit.



*Result:*



**Step**      **Action**

<b>UV monitor</b>	<b>Clamp placement</b>	<b>Result</b>
<b>U9-T</b> and <b>U9-L</b>	Place the base clamp to the left of the column.	 A photograph of a laboratory instrument's UV monitor assembly. A vertical white column is mounted on a base. A black clamp is attached to the left side of the column. At the top of the column, there is a red cap and a black sensor. The background shows a white laboratory bench with a 'Conductivity' label and a 'UV' label.
<b>U9-M</b>	Place the base clamp to the right of the column.	 A close-up photograph of the U9-M UV monitor assembly. A vertical white column is mounted on a base. A black clamp is attached to the right side of the column. The background shows a white laboratory bench with a 'U9-M' label.



Page intentionally left blank

**Give feedback on this document**

Visit [cytiva.com/techdocfeedback](https://cytiva.com/techdocfeedback) or scan the QR code.



## cytiva.com/akta

Cytiva and the Drop logo are trademarks of Life Sciences IP Holdings Corp. or an affiliate doing business as Cytiva.

ÄKTA, ÄKTA go, ÄKTA pure are trademarks of Global Life Sciences Solutions USA LLC or an affiliate doing business as Cytiva.

Any other third-party trademarks are the property of their respective owners.

© 2020–2022 Cytiva

For local office contact information, visit [cytiva.com/contact](https://cytiva.com/contact)

28968641 AG V:11 12/2022