

Research use only. Not for use in diagnostic procedures.

## Technical datasheet: Dimensions of DELFIA Streptavidin-coated 384-well microplate (white)

---

<b>Part numbers:</b>	PN CC11-H10	White, Streptavidin-coated
	PN CC11-H10S	White, Streptavidin-coated (Sample)
<b>Material and treatment:</b>	Polystyrene 384-well white microplate coated with streptavidin.	
<b>Packaging:</b>	CC11-H10: Contains 10 pcs of Streptavidin-coated 384-well plates in pouches. CC11-H10S: Sample pack of 2 pouched plates	
<b>Intended Use:</b>	The DELFIA Streptavidin-coated 384 well plates are intended for <i>in vitro</i> use only, as a solid phase in binding assays, e.g. fluoro-immunoassays.	
<b>Storage:</b>	Store plates at 2-8°C. NOTE: make sure the pouch remains sealed during storage to avoid moisture!	
<b>Coating Volume:</b>	The wells of this plate have been passively coated with 60µL of streptavidin and saturated with 80µL bovine serum albumin (BSA).	
<b>Warnings and Precautions:</b>	The DELFIA Streptavidin-coated 384 well plates are intended for research use only. Disposal of all waste should be in accordance with local regulations.	

### Warranty

Purchase of the product gives the purchaser the right to use this material in his own research, development, and investigational work. The product is not to be used for diagnostic procedures. Revvity, Inc. reserves the right to discontinue or refuse orders to any customer who plans to use products for any other purposes.

Revvity, Inc. does not warrant or guarantee that the product is merchantable or satisfactory for any particular purpose, nor free from any claim of foreign or domestic patent infringement by a third party, and there are no warranties, expressed or implied, to such effect.

Revvity, Inc. will not be liable for any incidental, consequential or contingent damages involving their use including damages to the property or personal injuries.

All information supplied with the product and technical assistance given is believed to be accurate, but it remains the responsibility of the investigator to confirm all technical aspects of the application. We appreciate receiving any additions, corrections, or updates to information supplied to the customer.

Dimensions -- Customer Drawing:

