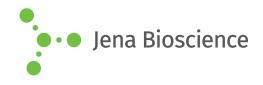
DATA SHEET





HDP Hanging Drop Crystallization Plate

material UVP

Cat. No.	Amount
CPL-151	10 pcs.

SWISSCI





For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: n/a

Form: box of 10 plates

The Swissci Hanging Drop Crystallization Plate

The Swissci Hanging Drop Crystallization Plate is a breakthrough for protein crystallization in a 96-well format. For the first time individual wells can be identified and removed without any disturbance to the growing crystals inside the plate.

The plate has been developed by Swissci AG together with leading Crystallographers. It is the result of many years of experience in successful robotic high-throughput crystallization.

The Hanging Drop Crystallization Plate offers unique properties that make it ideal for both nanoliter crystallization screening and microliter optimization alike. Made from optically superior polymer (UVP) and with a new design of the wells, the plate allows easy crystal viewing and retrieval.

The system is fully covered by design and patent protection.

Advantages of the Hanging Drop Crystallization Plate:

Easy crystal retrieval

Raised wide wells make the crystal mounting especially easy. Allen key system allows for smooth removal and subsequent microscopic investigation / X-ray with the novel crystal plug closure all included in the system.

Easy viewing

The wells are wide conical and have a lens effect for perfect illumination.

The micro-numbering ensures you will never loose the location under the microscope.

The optically superior polymer is UV transmissible and can be used to differentiate between salt and protein crystals.

Integral sealing

Wide partition walls between each well provide a high surface area for very good sealing with the advanced quality sealing tape. The sealing tape is integrated into the kit so there is no need to purchase additional parts prior to starting a procedure.

Wide range of volumes

Typical volumes are 50-200 µl of reservoir and 10 nl-5 µl drop size.

SBS standard

The complete plate is designed to the 96-well SBS standard for all common holders.

