



### MRC Maxi Plate - UVP

48-well plate  
highly UV-transparent polymer

Cat. No.	Amount
CPL-154S	10 pcs.
CPL-154L	100 pcs.




**For general laboratory use.**

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at ambient temperature

**Shelf Life:** n/a

**Form:** Plates wrapped inverted in sets of 10 double-bagged, clean room produced.

#### MRC Maxi 48-well Crystallisation Plate for Automated Optimisation

The MRC Maxi optimisation plate is a breakthrough for macromolecular crystallisation presented in a 48 well format. Offering easy to automate crystallisation optimisation with large sitting-drops, the new MRC Maxi Crystallisation plate is the perfect solution. Manufactured by Swissci AG, the plate offers an SBS format while providing 48 wells. MRC Maxi is intended for large drops and is compatible both with standard robotic systems as well as manual pipetting.

The plate was developed at the MRC Laboratory of Molecular Biology (Cambridge, UK) in collaboration with Jan Lowe and Fabrice Gorrec. It is a result of many years of experience in successful robotic high-throughput crystallisation and complements the original MRC crystallisation plate, which is intended for smaller drop volumes and higher throughput during screening.

Drop volumes of up to 10  $\mu$ l are possible. The 9 mm standard distance between wells is preserved, enabling the use of multi-channel manual pipettes and robotic liquid handlers, making MRC Maxi one of the most automation-friendly optimisation plates on the market.

MRC Maxi is covered by global intellectual property and design registration as are the Swissci AG MRC 2 lens 96 well plates. Several breakthrough features of the original MRC plate have been maintained. Wells are labeled individually. Drops are raised for easy access during crystal retrieval. MRC Maxi uses the same proprietary polymer specially selected for the purpose of UV light visualisation and the material used reduces through-plastic evaporation to a minimum. Well shapes are spherical but shallow.

#### The advantages of the MRC Maxi Crystallisation plate - in brief

##### Easy crystal retrieval

Raised wide wells make the crystal mounting especially easy.

##### Easy viewing

The wells are wide conical.

Each well has a micro lens for perfect illumination.

Micro numbering readable under the microscope for each well.

The optically superior polymer (UVP) is UV transmissible.

##### Better sealing

Wide partition walls between the wells give plenty of area for good sealing with tape.

Very rigid, automation-friendly plate design.

The UVP polymer reduces through-plastic evaporation to a minimum.

##### SBS Standard

The plates are designed to the SBS standard and are compatible with all common holders.

9 mm distance from well-to-well within columns, 18 mm distance within rows.



## MRC Maxi Plate - UVP

48-well plate  
highly UV-transparent polymer

Unique Polymer (UVP)  
Ultra-low sample binding.  
No static charging.

### Recommended volumes

Volumes validated for MRC Maxi are up to 10 µl of sample drop and 200 µl of the crystallisation reagent.