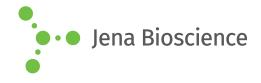
DATA SHEET

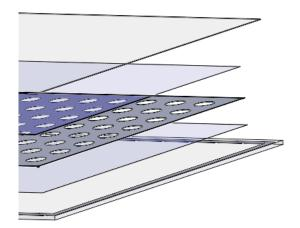




■ IMISX[™] Plate Kit

Includes components for 20 IMISX plates

Cat. No.	Amount
CPL-162	20 pcs.



IMISX™ Plate Layer View

For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: n/a

Description:

The IMISXTM Plate is designed to perform in meso in situ serial X-ray crystallography of soluble and membrane proteins [1].

The 96-well dual-sandwich IMISX™ Plate can be easily setup by hand or with an LCP dispensing robot. An SBS standard 96-well microplate footprint makes this plate compatible with standard robots and imagers.

By combining the standard LCP crystallization plate format with an advanced X-ray transparent thin-film design, the IMISX™ Plate allows for crystal growth and X-ray diffraction data collection like never before.

Content:

- 20 Glass LCP base plates with perimeter gaskets installed
- 20 Glass LCP top cover plates
- 40 Ultra Thin LCP Films (with protective liners) for inner sandwiches
- 20 96-well adhesive gaskets for LCP Film sandwiches
- 1 in situ Goniometer Adaptor & holder
- 5 Goniometer Adaptor mounting gaskets
- 1 Bottle of Silanization fluid
- 6 Clean wipe applicators
- Instructions

Technical Specifications

Base Plate (outer sandwich bottom)

Material: Glass Thickness: 1.1 mm Length: 127.7 mm Width: 85.5 mm

Notes: SBS standard, works with robots and storage hotels

Thin Film (inner sandwich top and bottom)
Material: UV and X-ray transparent Polymer

Thickness: 25 µm Length: 112 mm Width: 77 mm

Top Plate (outer sandwich top)

Material: Glass Thickness: 150 μm Length: 124 mm Width: 88 mm

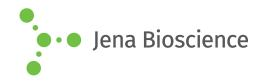
96-Well Adhesive Gasket

Material: Double-sided adhesive gasket

Well diameters: 5.0 mm Thickness: 140 µm Length: 112 mm Width: 77 mm



DATA SHEET





IMISX™ Plate Kit

Includes components for 20 IMISX plates

Selected References:

[1] Huang et al. (2015) In meso in situ serial X-ray crystallography of soluble and membrane proteins. Acta Cryst D 71:1238.

nonsaleable to USA and Japan