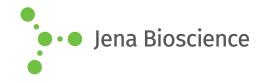
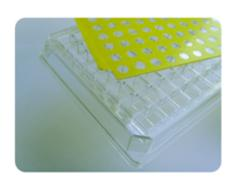
DATA SHEET





Lipidic Cubic Phase screening kit

Cat. No.	Amount
CPL-156	20 plates







For general laboratory use.

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: n/a

Form: box of 20 plates

Description:

The product comprises a base plate with low tack 700 micron thickness slide and top sealing tape of 100 microns - a brown ready to remove cover sheet and separate 200 micron UVP cover film with protective dust cover.

The LPC Kit

The Lipidic Cubic Phase Crystallizations (LCP) kit facilitates the automation and increased throughput of LCP crystallization set-ups. This novel system enables LCP screening to be performed accurately and with ease - using manual - or automated systems to complete the delivery of the solutions.

Working with innovative scientists Professor Gebhard Schertler and Pat Edwards from the MRC laboratory in Cambridge, Swissci AG developed a new system for setting up LCP screens.

The Swissci LCP kit ensures crystallography set-up is fast, cost effective and robotically adaptable. In addition Ultra Violet light can be used with the glass / UVP polymers to visualize proteins without the light scattering associated with traditional polymeric solutions.

The LCP technique for crystallising membrane proteins can be difficult and time-consuming to set up as it utilises highly viscous lipid mesophases to reconstitute proteins. Swissci AG and MRC overcame these problems with the LCP kit. This product combines an easy to use sandwich system with dedicated coupling tapes and thin film UVP cover plates. It conforms to SBS standards and thus is perfect for modern automation applications where high throughput is required.

The advantages of the LCP kit

- The plate design complies with SBS standard and prevents from cross contamination.
- Unique polymer: UV visualization is not compromised by polarization.
- Easy sealing with dry tabbed adhesive tape exposure and thin UVP cover film.

DATA SHEET





Lipidic Cubic Phase screening kit

 Unique low tack plate security allows for the sandwich plate to be removed from the base plate when required - in-situ X-ray data collection and structure determination is then enabled.