

**JBScreen Classic 1**

based on PEG 400 to 3000

Cat. No.	Amount
CS-101L	24 solutions (10 ml each)

For general laboratory use.**Shipping:** shipped at ambient temperature**Storage Conditions:** store at 8-10 °C**Shelf Life:** 12 months**Description:**

JBScreen Classic is a crystallization kit designed for efficient and flexible screening of crystallization conditions for proteins, peptides, nucleic acids, macromolecular complexes and water-soluble small molecules.

The JBScreen Classic Kits 1-10 cover 240 of the most prominent buffers for protein crystallization. Their compositions result from data mining of several thousands of crystallized proteins. JBScreen Classic represents the statistically most successful buffers that yielded protein crystals suitable for X-ray diffraction.

Content:

Set of 24 tubes in 10 ml bulk format with buffers based on PEG 400 to 3000.

The exact composition of each buffer is listed under the links "Screen Formulation".

All solutions are sterile filtered (0.2 µm), filled under sterile conditions and contain 0.1% ProClin™ 950 as preservative to prevent the growth of microbes. (ProClin is a trademark of LANXESS Corp.)

Related Products:

All solutions are available in customized formulations or larger quantities.