

**JBScreen Wizard 4**

Cat. No.	Amount
CS-314	48 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 Wizard is a highly effective random sparse matrix screen for crystallizing proteins, peptides, nucleic acids and macromolecular complexes. A large range of precipitants, buffers and salts allow a broad sampling of crystallization space at pH levels from pH 4,5 to 10,5.

**Content:**

Set of 48 tubes in 10 ml bulk format with unique screening solutions. 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.