

**JBScreen Plus 1**

kosmotropic reagents

Cat. No.	Amount
CS-501	24 solutions (1,0 ml each)

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

JBScreen Plus is an additive screen most useful in the optimization of preliminary crystallization conditions. The selection of the additives is based on the Hofmeister series, which reflects the ability of ions to stabilize the structure of proteins. Thus ions can be classified as either kosmotropic or chaotropic. The first having structure stabilizing properties, thus they may assist in, e.g. crystallizing proteins with a high proportion of flexible loop regions. The latter show structure disturbing properties which may assist in the crystallization of large complexes allowing them to re-arrange to form favorable crystal contacts.

Content:

Set of 24 tubes with 1 ml filling volume from a selection of kosmotropic reagents.

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