

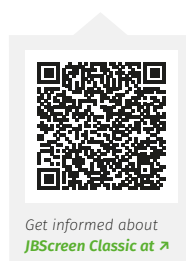
JBScreen Classic is a crystallization kit designed for efficient and flexible screening of crystallization conditions for proteins, peptides, nucleic acids, macromolecular complexes and water-soluble small molecules.

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.

The JBScreen Classic buffers are principally **ordered by type and concentration of the precipitant**. This allows easy extraction of all relevant information and is already a first step to a refinement: **Once you get a hit, you immediately see the effects of the neighbouring conditions**. Subsequent fine tuning of preliminary hits will be much more efficient.

JBScreen Classic comprises **10 kits of 24 unique reagents in the standard 10 ml bulk format**. JBScreen Classic HTS I+II contains the formulations of the JBScreen system, adopted to fit the 96-well format for high throughput crystallization applications. Each JBScreen Classic **HTS deep-well block is pre-filled with 96 sterile conditions at 1.7 ml each**.

Product	Cat-No.	Amount
JBScreen Classic 1 (PEG 400 to 3000 based)	CS-101L	24 solutions (10 ml each)
JBScreen Classic 2 (PEG 4000 based)	CS-102L	24 solutions (10 ml each)
JBScreen Classic 3 (PEG 4000+ based)	CS-103L	24 solutions (10 ml each)
JBScreen Classic 4 (PEG 5000 MME to 8000 based)	CS-104L	24 solutions (10 ml each)
JBScreen Classic 5 (PEG 8000 to 20000 based)	CS-105L	24 solutions (10 ml each)
JBScreen Classic 6 (Ammonium Sulfate based)	CS-106L	24 solutions (10 ml each)
JBScreen Classic 7 (MPD based)	CS-107L	24 solutions (10 ml each)
JBScreen Classic 8 (MPD/Alcohol based)	CS-108L	24 solutions (10 ml each)
JBScreen Classic 9 (Alcohol/Salt based)	CS-109L	24 solutions (10 ml each)
JBScreen Classic 10 (Salt based)	CS-110L	24 solutions (10 ml each)
JBScreen Classic 1 – 5	CS-112L	5 Kits
JBScreen Classic 6 – 10	CS-113L	5 Kits
JBScreen Classic 1 – 10	CS-114L	10 Kits
JBScreen Classic HTS I (PEG based)	CS-201L	96 solutions (1,7 ml each)
JBScreen Classic HTS II (Ammonium Sulfate, MPD, Alcohol and Salt based)	CS-202L	96 solutions (1,7 ml each)





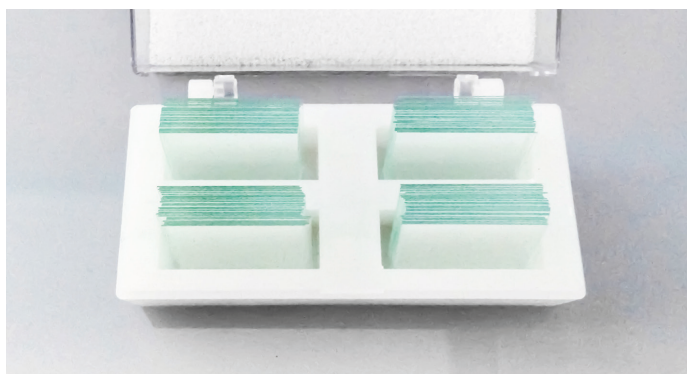
SUPERCLEAR 24 WELL PLATES

- › Made of crystal clear polystyrene for perfect visualization of droplets
- › Available pregreased or non-greased
- › Stackable, lid included
- › Individually wrapped
- › To be sealed with 22 mm cover slides



Get informed about
[SuperClear Plates](#) ↗

Product	Cat-No.	Amount
Crystalgen SuperClear Plates	CPL-130	100 pcs.
Crystalgen SuperClear Plates, pregreased	CPL-132	100 pcs.



COVER SLIDES

Glass cover slides are available plain or with a siliconized surface. 18 mm cover slides fit the Linbro Plate Space Saver, while 22 mm cover slides fit the SuperClear plates. Ideally suited for all kinds of crystallization experiments such as sandwich, hanging and sitting drop.



Get informed about
[Cover Slides](#) ↗

Product	Cat-No.	Amount
18 mm Circular Cover Slides – plain circle diameter: 18 mm; thickness: 0,22 mm	CSL-103	1000 pcs.
22 mm Circular Cover Slides – plain circle diameter: 22 mm; thickness: 0,22 mm	CSL-104	1000 pcs.
18 mm Circular Cover Slides – siliconized circle diameter: 18 mm; thickness: 0,22 mm	CSL-105	500 pcs.
22 mm Thick Circular Cover Slides – siliconized circle diameter: 22 mm; thickness: 0,5 mm	CSL-107	500 pcs.
18 mm Square Cover Slides – plain thickness: 0,22 mm	CSL-108	1000 pcs.

