

JBScreen Classic 3 (PEG 4000 + based)

The JBScreen Crystal Screening 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 represents the statistically most successful buffers that yielded protein crystals suitable for X-ray diffraction. The JBScreen 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. Apart from saving you the tedious pipetting, sterile packed aliquots have several other important advantages: While common bulk solutions are frequently contaminated by the time you use the first aliquot, JBScreen is always fresh and sterile. Evaporation of solvent (with the resulting change of the buffer) just won't happen - you will get extremely reproducible screening results.

The kit contains the following buffers:

- A1 8% PEG 4000, 0.8 M Li Chloride, 0.1 M TRIS HCl pH 8.5
- A2 10% PEG 4000, 20% iso-Propanol
- A3 10% PEG 4000, 10% iso-Propanol, 0.1 M Na Citrate pH 5.6
- A4 10% PEG 4000, 5% iso-Propanol, 0.1 M Na HEPES pH 7.5
- A5 10% PEG 4000, 20% iso-Propanol, 0.1 M Na HEPES pH 7.5
- A6 12% PEG 4000, 0.1 M Na Acetate pH 4.6, 0.2 M Ammonium Sulfate
- B1 15% PEG 4000, 0.2 M Ammonium Sulfate
- B2 15% PEG 4000, 0.1 M Na Citrate pH 5.6, 0.2 M Ammonium Sulfate
- B3 16% PEG 4000, 10% iso-Propanol, 0.1 M Na HEPES pH 7.5, 0.2 M Ammonium Sulfate
- B4 20% PEG 4000, 0.2 M Ammonium Sulfate
- B5 20% PEG 4000, 10% Glycerol, 0.2 M Mg Sulfate
- B6 20% PEG 4000, 5% iso-Propanol, 0.1 M Na Citrate
- C1 20% PEG 4000, 20% iso-Propanol, 0.1 M Na Citrate
- C2 20% PEG 4000, 0.6 M Na Chloride, 0.1 M Na MES pH 6.5
- C3 20% PEG 4000, 10% iso-Propanol, 0.1 M Na HEPES pH 7.5
- C4 22% PEG 4000, 0.2 M Ammonium Sulfate, 0.1 M Na Acetate
- C5 25% PEG 4000, 0.1 M Na Acetate pH 4.6, 0.2 M Ammonium Sulfate
- C6 25% PEG 4000, 0.1 M Na Citrate pH 5.6, 0.2 M Ammonium Sulfate
- D1 25% PEG 4000, 0.1 M Na HEPES pH 7.5, 0.2 M Li Sulfate, 0.1 M Na Acetate

D2 25% PEG 4000, 8% iso-Propanol, 0.1 M Na Acetate

D3 30% PEG 4000, 0.2 M Ammonium Sulfate

D4 30% PEG 4000, 0.1 M Na Acetate pH 4.6, 0.2 M Ammonium Sulfate

D5 30% PEG 4000, 0.1 M Na Citrate pH 5.6, 0.2 M Ammonium Sulfate

D6 32% PEG 4000, 0.8 M Li Chloride 0.1 M TRIS HCl pH 8.5

Buffer preparation:

- Chemicals used were of MicroSelect grade for Molecular Biology.
- Buffers were prepared as 1 M stock solutions. The pH was adjusted with HCl (Imidazole, TRIS), NaOH (Bicine, CHES, HEPES, MES), Citric Acid (Citrate) or Acetic Acid (Acetate) (23°C; Fisher pH electrode).
- Percentages given are always w/v values (for solids as well as for liquids).
- The final volume was adjusted with >18 MOhm water.
- Solutions were filtered (0.2 µm) and filled in ampoules/tubes under sterile conditions.

Ordering Information

[Order Online »](#)

JBS-CS-103L	1Kit
-------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Contents

JBS-CS-103: 24 buffers, delivered in 0.7ml sterile aliquots, with 10 units per kit. Alternatively, the JBScreen Mixed kit contains one single shot of all the 240 buffers. The 6×4 format of the units allows a convenient 1:1 setup for 24 (or multiples thereof) well plates. It's easy to use – every ampoule and the unit's package has a “coordinate system” (A1–A6 ... D1–D6), thus almost certainly eliminating errors.

JBS-CS-103L: All 24 solution of the JBScreen system in sterile 10ml quantities (Bulk).

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)