

# JBScreen Classic 1

## (PEG 400 to 3000 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 15% PEG 400, 0.1 M Na Acetate pH 4.6, 0.1 M Ca Chloride
- A2 15% PEG 400, 0.1 M Na MES pH 6.5
- A3 15% PEG 400, 0.1 M Na HEPES pH 7.5, 0.2 M Mg Chloride
- A4 15% PEG 400, 0.1 M TRIS HCl pH 8.5, 0.2 M Na Citrate
- A5 25% PEG 400, 0.1 M Na Acetate pH 4.6, 0.1 M Mg Chloride
- A6 25% PEG 400, 0.1 M TRIS HCl pH 8.5, 0.2 M Li Sulfate
- B1 28% PEG 400, 0.1 M Na HEPES pH 7.5, 0.2 M Ca Chloride
- B2 30% PEG 400, 0.1 M Na Acetate pH 4.6, 0.1 M Ca Chloride
- B3 30% PEG 400, 0.1 M Na MES pH 6.5, 0.1 M Na Acetate
- B4 30% PEG 400, 0.1 M Na MES pH 6.5, 0.1 M Mg Chloride
- B5 30% PEG 400, 0.1 M Na HEPES pH 7.5, 0.2 M Mg Chloride
- B6 30% PEG 400, 0.1 M TRIS HCl pH 8.5, 0.2 M Na Citrate
- C1 30% PEG 550 MME, 0.1 M Bicine pH 9.0, 0.1 M Na Chloride
- C2 25% PEG 550 MME, 0.1 M Na MES pH 6.5, 0.01 M Zn Sulfate
- C3 25% PEG 1000, 0.1 M Na HEPES pH 7.5
- C4 30% PEG 1000, 0.1 M TRIS HCl pH 8.5
- C5 15% PEG 1500
- C6 20% PEG 1500, 0.1 M Na HEPES pH 7.5
- D1 30% PEG 1500
- D2 20% PEG 2000 MME, 0.1 M TRIS HCl pH 8.5, 0.01 M Ni Chloride
- D3 25% PEG 2000 MME

D4 30% PEG 2000 MME, 0.1 M Na MES pH 6.5, 0.1 M Na Acetate

D5 20% PEG 3000, 0.1 M Na HEPES pH 7.5, 0.2 M Na Acetate

D6 30% PEG 3000, 0.1 M TRIS HCl pH 8.5, 0.2 M Li Sulfate

## Ordering Information

[Order Online »](#)

JBS-CS-101L	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-101:** 24 buffers, delivered in 0.7 ml 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-101L:** All 24 solution of the JBScreen system in sterile 10ml quantities (Bulk).

**JBS-CS-101S:** 24×10 sterile aliquots of the JBScreen buffers, supplied as 10-ampoule blocks, with 0.7ml per ampoule (Refill).

**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)