

# JBScreen Plus

## Kosmotropic

JBScreen Plus solutions are most useful in the optimization of preliminary crystallization conditions, where specific interactions between the additives and the biological macromolecule may assist in the initial crystallization of the sample and may help to improve the size and diffraction quality of the crystals.

JBScreen Plus comprises 5 kits, including kosmotropic (structure-stabilizing) and chaotropic (structure-disturbing) additives, salts, volatile and non-volatile organics and other compounds. Each kit contains 24 different additives, supplied in ready-made aliquots (100 µl each, except the JBScreen Plus Volatiles kit, containing 1 ml each), with adjusted concentration and sterile filtered. All solutions come in tightly closing glass vials that will keep them fresh and sterile.

### Ordering Information

[Order Online »](#)

JBS-CS-501	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 24 different compounds, 100µl each, delivered in 2ml screw top vials.

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)