

JBScreen Plus

Additives

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-504	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 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

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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