

JBScreen Basic 3

In 1991, Jancarik and Kim published 50 conditions, which were derived from previously crystallized proteins [1]. These and other conditions form the basis of the JBScreen Basic system. However, JBScreen Basic is designed to fit the 24-well plate format and like in all other JBScreen crystallization kits, we abstained from the use of cacodylate buffers and replaced them with MES. JBScreen Basic contains 96 unique reagent mixtures for screening a wide range of pH and various salts and precipitants. Each condition of the four kits is supplied in 10 ml quantities.

Ordering Information

[Order Online »](#)

JBS-CS-123	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 unique, sterile filtered reagents, supplied in 10ml aliquots.

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