

JBScreen Classic HTS II (Ammonium Sulfate, MPD, Alcohol and Salt based)

JBScreen HTS is designed for high throughput crystallization applications using multi-channel pipets or crystallization robots. Each JBScreen HTS 96-well masterblock is pre-filled with 96 sterile crystallization buffers, with either 1.0 ml (S size) or 1.7 ml (L size) per well and is sealed with indexed cap mats.

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.

For a detailed list of buffers contained in the kit please [click here](#):

Ordering Information

[Order Online »](#)

JBS-CS-202L	1Kit
-------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

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