

JBScreen Membrane 3 (Ammonium Sulfate/Alcohol/Salt based)

The JBScreen Membrane Screens 1-3 cover 72 of the most successful buffers for crystallization of membrane proteins. Their compositions result from analyzing the crystallization conditions of all membrane proteins crystallized so far.

The JBScreen Membrane crystallization buffers are principally ordered by type and concentration of the precipitant. Like in case of the "classic" JBScreen system, 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.

The kit contains the following buffers:

- A1 0.7 M Ammonium Sulfate, 1 M Na/K Phosphate pH 7.5
- A2 1 M Ammonium Sulfate, 1 % Heptane-1,2,3-triol, 100 mM Zn Acetate, 50 mM Na MES pH 6.5
- A3 1.2 M Ammonium Sulfate, 3 % Heptane-1,2,3-triol, 50 mM Tris HCl pH 7.5
- A4 1.2 M Ammonium Sulfate, 100 mM Tris HCl pH 8.5
- A5 1.4 M Ammonium Sulfate, 4 % iso-Propanol, 100 mM Ammonium Acetate
- A6 2 M Ammonium Sulfate
- B1 2 M Ammonium Sulfate, 100 mM Na Citrate pH 5.6
- B2 2.5 M Ammonium Sulfate, 2 % PEG 5000 MME, 100 mM Na HEPES pH 7.5
- B3 3 M Ammonium Sulfate, 3 % Heptane-1,2,3-triol
- B4 3.5 M Ammonium Sulfate
- B5 3.5 M Ammonium Sulfate, 250 mM Na Chloride, 50 mM Na/K Phosphate pH 7.5
- B6 25 % MPD, 100 mM Bis-Tris pH 7.0
- C1 30 % MPD, 300 mM Na Citrate pH 5.6
- C2 25 % Triethylene Glycol, 100 mM Ammonium Sulfate, 100 mM Glycine
- C3 30 % iso-Propanol, 20 % Glycerol, 200 mM Ca Chloride, 100 mM Na Acetate pH 4.6
- C4 50 mM K Phosphate pH 8.0
- C5 100 mM Na Citrate pH 4.8
- C6 1 M K Phosphate pH 6.5, 3 % Heptane-1,2,3-triol, 1 % 1,4-Dioxane
- D1 1 M Na Citrate

D2 1 M Na Citrate pH 5.6, 500 mM Li Chloride

D3 1.5 M Na Chloride, 100 mM Na Acetate pH 4.6

D4 1.5 M K Phosphate pH 7.0

D5 1.5 M Li Sulfate, 100 mM Na HEPES pH 7.5

D6 2 M Na Chloride, 100 mM Na Formate

Ordering Information

[Order Online »](#)

JBS-CS-303L	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 sterile solutions, 10ml each.

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