



Jena Bioscience, with over 25 years of experience, is a leading provider of inovative and high-quality reagents and customized services in the life science field.



Bulk – Tailored to your scale.

Ready-made or customized reagents in large quantities, with personalized packaging and scientific support.



OEM – Your products, your way.

Exclusivity with customized formulations, packaging and branding under strict confidentiality.

Contact our Xtals Experts:

Phone: +49(0)3641 - 6285 000
Mail: xtals@jenabioscience.com



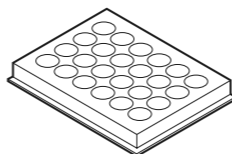
Dr. Bürk Schäfer
Head of Crystallography & Cryo-EM



PLATES & SEALINGS

24 Well Plates

- Linbro Plate
- SuperClear™ Plates
- ComboPlates™

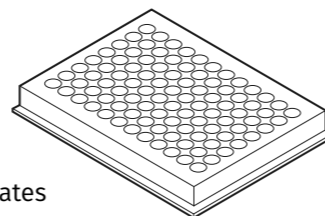


48 Well Plates

- MRC Maxi Plate

96 Well Plates

- In-situ-1 Plate
- CrystalDirect™ Plate
- MRC 2-well and 3-well Plates
- HDP Hanging Drop Plate



Lipidic Cubic Phase Plates

- IMISX™ Plate
- LCP Glass Sandwich Set
- Lipidic Cubic Phase Screening Kit
- HDP Hanging Drop Plate

Microbatch Plates

- MRC Under Oil Plate

Sealings

- Cover Slides & Grease, Sealing Tape

Miscellaneous

- 96 Well Micro-Dialysis Plate
- 96 Well Masterblocks & Cap Mats

CRYSTAL HANDLING

XtalTool

- Sample Holder for Crystal Growth, *in situ* Ligand Soaking and Data Collection

Retrieval and Mounting

- MicroMounts™, MicroLoops™, MicroMeshes™ and MicroGrippers™ from MiTeGen

Manipulation and Measurement

- MicroTool™ Kits from MiTeGen

Cryo and Room Temperature Crystallography

- Goniometer Bases (Caps), Magnetic CryoVials, Mount-Base-Vial Assemblies, Cryo Tools, MicroRT™ Tubing and MicroRT™ Aligner, Oils and Grease, Crystallography Starter Kits

Storage and Shipping

- Magnetic CryoVials, Pucks, Dry Shippers, Dewars and more...

Accessories

- Base Holders, Goniometer Head Adapter, Paper Wicks and Tweezers

YOUR PARTNER IN PROTEIN CRYSTALLIZATION



Screens
Optimization & Phasing
Plates & Sealings
Crystal Handling



CRYSTAL SCREEN SERIES

JBScreen Classic

- designed for efficient and flexible screening of crystallization conditions

JBScreen Basic

- for screening a wide range of pH and various salts and precipitants

JBScreen JCSG++

- formulated by researchers from the Joint Center for Structural Genomics (JCSG)

JBScreen Kinase

- optimized for highly effective crystallization of kinases

JBScreen Nuc-Pro

- for preliminary crystallization conditions of nucleic acids and protein-nucleic acid complexes

JBScreen PACT++

- systematic pH, anion- and cation testing in the presence of PEG

JBScreen PEG/Salt

- enables the screening of PEG depending on ionic strength, ion type and pH

JBScreen Wizard

- allow a broad sampling of crystallization space at pH levels from pH 4,5 to 10,5

SPECIAL SCREENS

XP Screens

- Anderson-Evans polyoxotungstate promotes crystallization of most challenging targets and improves diffraction quality

ThermoFluor Screens

- Thermal Shift Assay for protein stability

Additive Screens

- selection of the additives based on the Hofmeister series

Buffer Screens

- standard buffers and buffers with extreme pH range

FORMOScreen®

- 2bind Antibody Formulation Screen

JBScreens are available as:

Bulk format

- 24 × 10 ml in single tubes (or set of 4 × 24 tubes)

HTS format

- 96 × 1,7 ml in a deep-well block

OPTIMIZATION

Chemical Environment

- Kits to increase protein solubility and stability

Thermodynamics / Kinetics

- Improve diffraction by crystal annealing, varying the solvent content or use your initial crystals for seeding

Membrane Proteins

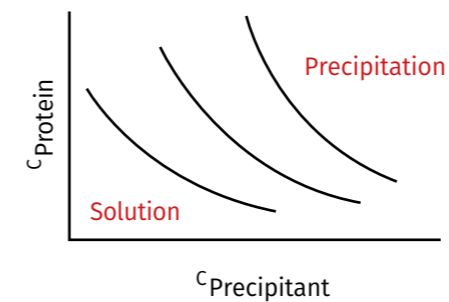
- LCP lipids & phospholipids to form a stable lipidic cubic phase for membrane protein crystallization

Crystal Dyes

- Choose the dye to stain your crystals blue, purple, green or red

Model Proteins

- Lysozyme solution as a model protein in crystallization quality



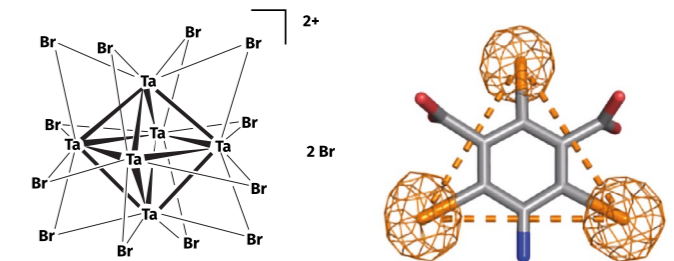
PHASING

Tantalum Cluster Derivatization Kit

JBS Tungsten Cluster Derivatization Kits

JBS Magic Triangle

Mercurated or Selenium-containing Nucleotides



Tantalum Bromide Cluster

JBS Magic Triangle

Carefully selected kits containing heavy atoms suitable for experiments involving single/multiple isomorph replacement (SIR and MIR) and anomalous dispersion (SAD and MAD).