

# JBScreen Classic – Crystallization Screening of Biological Macromolecules

**An efficient and flexible crystallization screening system for proteins, peptides, nucleic acids, macromolecular complexes and water-soluble small molecules.**



- **The JBScreen Classic crystallization screening kits represent a comprehensive statistical compilation of the most successful crystallization conditions from the thousands published in the Biological Macromolecule Crystallization Database [1].**
- **The full JBScreen Classic comprises 10 screening kits covering 240 conditions, with each kit containing reagent formulations grouped according to precipitant type and concentration.**
- **A broad sampling of crystallization space and select formulations proven to maximize the rate of success make JBScreen Classic the number one choice for academic and industrial labs.**

[1] Gilliland *et al.* (2002) The Biological Macromolecule Crystallization Database: crystallization procedures and strategies. *Acta Cryst. D* **58**:916.

## Our high quality crystal screens...



**JBScreen Classic bulk**



**JBScreen Classic HTS**

<b>Kit</b>	<b>Primary Precipitant</b>
JBScreen Classic 1	PEG 400 to 3000
JBScreen Classic 2	PEG 4000
JBScreen Classic 3	PEG 4000 +
JBScreen Classic 4	PEG 5000 MME to 8000
JBScreen Classic 5	PEG 8000 to 20000
JBScreen Classic 6	Ammonium Sulfate
JBScreen Classic 7	MPD
JBScreen Classic 8	MPD / Alcohol
JBScreen Classic 9	Alcohol / Salt
JBScreen Classic 10	Salt

## ...bring out the best in your project.

- **The largest, most comprehensive crystallization screen on the market**
- **Rational zone-organized layout of crystallization conditions**
- **Potent formulations based on decades of empirical data**
- **Two formats to satisfy all requirements**
  - **bulk: 10 individual kits @ 24x10 ml volumes**
  - **HTS: 2x96-well masterblocks for high-throughput applications (1.0 or 1.7 ml per well)**
- **Check our promotions for unbeatable prices!**

**Please contact us at [xtals@jenabioscience.com](mailto:xtals@jenabioscience.com)**

# Numerous literature citations show that JBScreen has been successfully employed in crystallization experiments

- Marcia *et al.* (2009) The structure of *Aquifex aeolicus* sulfide: quinone oxidoreductase, a basis to understand sulfide detoxification and respiration. *Proc. Natl. Acad. Sci.* **106**:9625.
- Dunstan *et al.* (2009) Structure of the Thiostrepton Resistance Methyltransferase\_S-Adenosyl-L-methionine Complex and Its Interaction with Ribosomal RNA. *J. Biol. Chem.*, **284**:17013.
- Sugishima *et al.* (2009) Crystal Structure of Red Chlorophyll Catabolite Reductase: Enlargement of the Ferredoxin-Dependent Bilin Reductase Family. *J. Mol. Biol.* **389**:376.
- Raaf *et al.* (2009) First Inactive Conformation of CK2 $\alpha$ , the catalytic Subunit of Protein Kinase CK2. *J. Mol. Biol.* **389**:1212.
- Cossu *et al.* (2009) Designing Smac-mimetics as antagonists of XIAP, cIAP1, and cIAP2. *Biochem. Biophys. Res. Commun.* **378**:162.
- Gourlay *et al.* (2009) Group B Streptococcus Pullulanase Crystal Structures in the Context of a Novel Strategy for Vaccine Development. *J. Bacteriology* **191**:3544.
- Jansson *et al.* (2009) Structure of the methyltransferase domain from the Modoc virus, a flavivirus with no known vector. *Acta Cryst. D* **65**:796.
- Souza *et al.* (2009) Crystallization, data collection and data processing of maltose-binding protein (MalE) from the phytopathogen *Xanthomonas axonopodis* pv. *citri*. *Acta Cryst. F* **65**:105.
- Cui *et al.* (2009) Crystallization and preliminary crystallographic analysis of the second RRM of Pub1 from *Saccharomyces cerevisiae*. *Acta Cryst. F* **65**:108.
- Prudnikova *et al.* (2009) Crystallization and preliminary X-ray analysis of a novel haloalkane dehalogenase DbeA from *Bradyrhizobium elkani* USDA94. *Acta Cryst. F* **65**:353.
- Tulsidas *et al.* (2009) Crystallization of a nonclassical Kazal-type *Carcinoscorpius rotundicauda* serine protease inhibitor, CrSPI-1, complexed with subtilisin. *Acta Cryst. F* **65**:533.
- Bogdanovic *et al.* (2009) Crystallization and preliminary X-ray diffraction studies of AsaP1\_E294A and AsaP1\_E294Q, two inactive mutants of the toxic zinc metallopeptidase AsaP1 from *Aeromonas salmonicida* subsp. *achromogenes*. *Acta Cryst. F* **65**:695.
- Squina *et al.* (2009) Expression, purification, crystallization and preliminary crystallographic analysis of an endo-1,5-L-arabinanase from hyperthermophilic *Thermotoga petrophila*. *Acta Cryst. F* **65**:902.
- Kudlinzki *et al.* (2009) Crystallization and preliminary X-ray diffraction analysis of the C-terminal domain of the human spliceosomal DExD/H-box protein hPrp22. *Acta Cryst. F* **65**:956.

- Weber *et al.* (2009) Crystallization and preliminary X-ray analysis of a D-Ala:D-Ser ligase associated with VanG-type vancomycin resistance. *Acta Cryst. F* **65**:1024.
- Reiser *et al.* (2009) Crystallization and preliminary X-ray crystallographic characterization of a public CMV-specific TCR in complex with its cognate antigen. *Acta Cryst. F* **65**:1157.
- Piton *et al.* (2009) Purification, crystallization and preliminary X-ray diffraction experiments on the breakage-reunion domain of the DNA gyrase from *Mycobacterium tuberculosis*. *Acta Cryst. F* **65**:1182.
- Vulliez-Le Normand *et al.* (2008) Structures of synthetic O-antigen fragments from serotype 2a *Shigella flexneri* in complex with a protective monoclonal antibody. *Proc. Natl. Acad. Sci.* **105**:9976.
- Kurz *et al.* (2008) Crystallization and preliminary diffraction analysis of a DsbA homologue from *Wolbachia pipientis*. *Acta Cryst. F* **64**: 94.
- Stsiapanava *et al.* (2008) Crystals of DhaA mutants from *Rhodococcus rhodochrous* NCIMB 13064 diffracted to ultrahigh resolution: crystallization and preliminary diffraction analysis. *Acta Cryst. F* **64**:137.
- Mueller-Dieckmann *et al.* (2008) Structure of mouse ADP-ribosylhydrolase 3 (mARH3). *Acta Cryst. F* **64**:156.
- Yamada *et al.* (2008) Crystallization and preliminary crystallographic analysis of the transpeptidase domain of penicillin-binding protein 2B from *Streptococcus pneumoniae*. *Acta Cryst. F* **64**:284.
- Pérez de Diego *et al.* (2008) Crystallization and preliminary X-ray diffraction analysis of the subunit Yke2 of the Gim complex from *Saccharomyces cerevisiae*. *Acta Cryst. F* **64**:501.
- Okada *et al.* (2007) Crystal Structure of the  $\gamma$ -Glutamyltranspeptidase Precursor Protein from *Escherichia coli*. *J. Biol Chem.* **282**:2433.
- Sato *et al.* (2007) Crystal structure of the Sec4p-Sec2p complex in the nucleotide exchanging intermediate state. *Proc. Natl. Acad. Sci.* **104**:8305.
- Cowan-Jacob *et al.* (2007) Structural biology contributions to the discovery of drugs to treat chronic myelogenous leukaemia. *Acta Cryst. D* **63**:80.
- Ingvarsson *et al.* (2007) Insights into the inter-ring plasticity of caseinolytic proteases from the X-ray structure of *Mycobacterium tuberculosis* ClpP1. *Acta Cryst. D* **63**:249.
- Hagemeyer *et al.* (2006) Insight into the mechanism of biological methanol activation based on the crystal structure of the methanol-cobalamin methyltransferase complex. *Proc. Natl. Acad. Sci.* **103**:18917.
- Smatanova' *et al.* (2006) New techniques for membrane protein crystallization tested on photosystem II core complex of *Pisum sativum*. *Photosynth Res* **90**:255.
- Dauter (2006) Current state and prospects of macromolecular Crystallography. *Acta Cryst. D* **62**:1.
- Rypniewski *et al.* (2006) The first crystal structure of an RNA racemate. *Acta Cryst. D* **62**:659.
- le Maire *et al.* (2006) Crystallization and halide phasing of the C-terminal domain of human KIN17. *Acta Cryst. F* **62**:245.
- Goyal *et al.* (2006) Crystallization and preliminary X-ray crystallographic studies of the N-terminal domain of FadD28, a fatty-acyl AMP ligase from *Mycobacterium tuberculosis*. *Acta Cryst. F* **62**:350.
- Tsukazaki *et al.* (2006) Purification, crystallization and preliminary X-ray diffraction of SecDF, a translocon-associated membrane protein, from *Thermus thermophilus*. *Acta Cryst. F* **62**:376.
- Ochiai *et al.* (2006) Crystallization and preliminary X-ray analysis of the rhamnogalacturonan lyase YesW from *Bacillus subtilis* strain 168, a member of polysaccharide lyase family 11. *Acta Cryst. F* **62**:438.

- Hilario *et al.* (2006) Crystallization and preliminary X-ray diffraction analysis of XAC1151, a small heat-shock protein from *Xanthomonas axonopodis* pv. *citri* belonging to the  $\alpha$ -crystallin family. *Acta Cryst. F* **62**:446.
- Matera *et al.* (2006) Preliminary crystallographic analysis of salicylate 1,2-dioxygenase from *Pseudaminobacter salicylatoxidans*. *Acta Cryst. F* **62**:553.
- Chuenchor *et al.* (2006) Purification, crystallization and preliminary X-ray analysis of rice BGlu1 b-glucosidase with and without 2-deoxy-2-fluoro- $\beta$ -D-glucoside. *Acta Cryst. F* **62**:798.
- Mastrangelo *et al.* (2006) Preliminary crystallographic characterization of an RNA helicase from Kunjin virus. *Acta Cryst. F* **62**:876.
- Stout *et al.* (2006) Crystallization, preliminary crystallographic analysis and phasing of the thiosulfate-binding protein SoxY from *Chlorobium limicola* f. *thiosulfatophilum*. *Acta Cryst. F* **62**:1093.
- Koch *et al.* (2006) Purification and crystallization of the human EF-hand tumour suppressor protein S100A2. *Acta Cryst. F* **62**:1120.
- Faravelli *et al.* (2006) Expression, refolding and crystallizations of the Grb2-like (GADS) C-terminal SH3 domain complexed with a SLP-76 motif peptide. *Acta Cryst. F* **62**:1278.
- Hülsmeyer *et al.* (2005) A major histocompatibility complex-peptide-restricted antibody and T cell receptor molecules recognize their target by distinct binding modes. *J. Biol. Chem.* **280**:2972.
- Cherezov *et al.* (2005) A simple and inexpensive nanoliter-volume dispenser for highly viscous materials used in membrane protein crystallization. *J. Appl. Cryst.* **38**:398.
- Judge *et al.* (2005) An ultraviolet fluorescence-based method for identifying and distinguishing protein crystals. *Acta Cryst. D* **61**:60.
- Miyatake *et al.* (2005) Development of a fully automated macromolecular crystallization/observation robotic system, HTS-80. *Acta Cryst. D* **61**:658.
- Leulliot *et al.* (2005) High-throughput crystal-optimization strategies in the South Paris Yeast Structural Genomics Project: one size fits all? *Acta Cryst. D* **61**:664.
- Huynh *et al.* (2005) Expression, purification and preliminary crystallographic analysis of sucrose phosphate synthase (SPS) from *Halothermothrix orenii*. *Acta Cryst. F* **61**:116.
- Mueller *et al.* (2005) Crystallization and preliminary X-ray diffraction analysis of human growth and differentiation factor 5 (GDF-5). *Acta Cryst. F* **61**:134.
- Park *et al.* (2005) Crystallization and preliminary X-ray crystallographic analysis of peptide deformylase (PDF) from *Bacillus cereus* in ligand-free and actinonin bound forms. *Acta Cryst. F* **61**:150.
- Ferraroni *et al.* (2005) Crystallization and preliminary structure analysis of the blue laccase from the ligninolytic fungus *Panus tigrinus*. *Acta Cryst. F* **61**:205.
- Schmidberger *et al.*, (2005) Purification, crystallization and preliminary crystallographic analysis of DehIVa, a dehalogenase from *Burkholderia cepacia* MBA4. *Acta Cryst. F* **61**: 271
- Ashikawa *et al.* (2005) Crystallization and preliminary X-ray diffraction analysis of the electron-transfer complex between the terminal oxygenase component and ferredoxin in the Rieske non-haem iron oxygenase system carbazole 1,9a-dioxygenase. *Acta Cryst. F* **61** :577.

- Harada *et al.* (2005) Crystallization and preliminary X-ray diffraction study of BchU, a methyltransferase from *Chlorobium tepidum* involved in bacteriochlorophyll *c* biosynthesis. *Acta Cryst. F* **61**:712.
- Lee *et al.* (2005) Crystallization and preliminary X-ray crystallographic analysis of human RGS10 complexed with Gai3. *Acta Cryst. F* **61**:831
- Beaven *et al.* (2005) Crystallization and preliminary X-ray diffraction analysis of calyculin from *Loligo pealei*: a neuronal protein implicated in learning and memory. *Acta Cryst. F* **61**:879.
- Songsirithigul *et al.* (2005) Expression, purification, crystallization and preliminary crystallographic analysis of chitinase A from *Vibrio carchariae*. *Acta Cryst. F* **61**:895.
- Irimia *et al.* (2004) Methanoarchaeal sulfolactate dehydrogenase: prototype of a new family of NADH-dependent enzymes. *EMBO J.* **23**:1234.
- Kundrot *et al.* (2004) Which strategy for a protein crystallization project? *Cell Mol. Life Sci.* **61**:525.
- Vallazza *et al.* (2004) First look at RNA in L-configuration *Acta Cryst. D* **60**:1.
- Unciuleac *et al.* (2004) Crystallization of 4-hydroxybenzoyl-CoA reductase and the structure of its electron donor ferredoxin. *Acta Cryst. D* **60**:388
- Sjekloca *et al.* (2004) Cloning, expression, purification, crystallization and preliminary crystallographic analysis of  $\gamma$ -filamin repeat 23. *Acta Cryst. D* **60**:1155.
- Cherezov *et al.* (2004) A robotic system for crystallizing membrane and soluble proteins in lipidic mesophases. *Acta Cryst. D* **60**:1795.
- Heras *et al.* (2003) Dehydration converts DsbG crystal diffraction from low to high resolution. *Structure* **11**:139.
- Vorup-Jensen *et al.* (2003) Structure and allosteric regulation of the  $\alpha\chi\beta 2$  integrin I domain. *Proc. Natl. Acad. Sci. USA* **100**:1873.
- Kimber *et al.* (2003) Data mining crystallization databases: Knowledge-based approaches to optimize protein crystal screens. *Proteins* **51**:562
- Smatanova *et al.* (2003) Proteins and their crystals. *Materials Structure* **10**:30.
- Georgieva *et al.* (2003) Crystallization and preliminary X-ray diffraction studies of a toxic phospholipase A2 from the venom of *Vipera ammodytes meridionalis* complexed to a synthetic inhibitor. *Biochim. Biophys. Acta* **1650**:1.
- Choi *et al.* (2003) Crystallization and preliminary X-ray analysis of *Escherichia coli* p20, a novel thiol peroxidase. *Acta Cryst. D* **59**:1064.
- Fermani *et al.* (2003) Crystallization and preliminary X-ray diffraction analysis of two ribosome-inactivating proteins: lychnin and dianthin 30. *Acta Cryst. D* **59**:1227.
- Kriksunov *et al.* (2003) Crystallization and preliminary crystallographic analysis of a new class of glutathione transferase from nematodes. *Acta Cryst. D* **59**:1262.
- Hagemeyer *et al.* (2003) Coenzyme F<sub>420</sub>-dependent methylenetetrahydromethanopterin dehydrogenase from *Methanopyrus kandleri*: the selenomethionine-labelled and non-labelled enzyme crystallized in two different forms. *Acta Cryst. D* **59**:1653.
- Makde *et al.* (2003) Expression, purification, crystallization and preliminary X-ray diffraction studies of recombinant class B non-specific acid phosphatase of *Salmonella typhimurium*. *Acta Cryst. D* **59**:1849.

- Dimasi *et al.* (2003) Expression, crystallization and preliminary crystallographic analysis of the extracellular IgV-like domain of the human natural killer cell inhibitory receptor p75/AIRM1. *Acta Cryst. D* **59**:1856.
- Kim *et al.* (2003). Crystallization and preliminary X-ray analysis of the small component of 4-hydroxyphenylacetate 3-monooxygenase (HpaC) and its cofactor complex from *Thermus thermophilus* HB8. *Acta Cryst. D* **59**, 2275.
- Fieulaine *et al.* (2002) X-ray structure of a bifunctional protein kinase in complex with its protein substrate HPr. *Proc. Natl. Acad. Sci. USA* **99**:13437.
- Larivière *et al.* (2002) Crystallization and preliminary crystallographic study of a ternary complex between the T4 phage  $\beta$ -glucosyltransferase, uridine diphosphoglucose and a DNA fragment containing an abasic site. *Acta Cryst. D* **58**:1484.
- Vallazza *et al.* (2002) Crystallization and structure analysis of *Thermus flavus* 5S rRNA helix B. *Acta Cryst. D* **58**:1700.
- Pesce *et al.* (2002) Human neuroglobin: crystals and preliminary X-ray diffraction analysis. *Acta Cryst. D* **58**:1848.