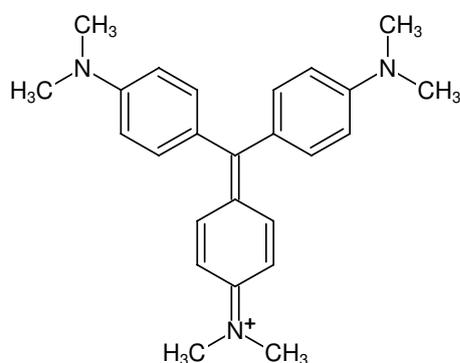


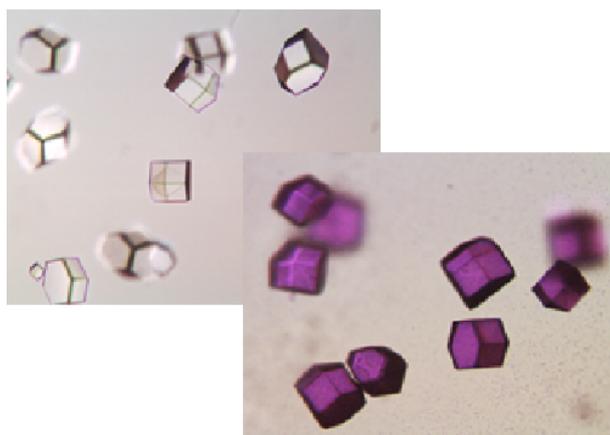

JBS Deep Purple

Crystal Violet

Cat. No.	Amount
CO-302	300 µl

Cl⁻

Structural formula of JBS Deep Purple



Unstained (A) and stained (B) protein crystals

For general laboratory use.
Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: 12 months

Molecular Formula: C₂₅H₃₀N₃Cl

Molecular Weight: 407.98 g/mol

CAS#: 548-62-9

EC number: 208-953-6

Concentration: 0.5 mM

Applications:

JBS Deep Purple is a crystal dye used to stain macromolecular crystals, i.e. protein, peptide and nucleic acid crystals in order to differentiate them from small molecules and salt crystals.

Description:

Crystallization screening with high concentrations of precipitant and salt may lead to the formation of salt crystals. It is quite difficult to make a distinction between these false positives and true protein crystals.

Staining of crystals with appropriate dyes is a very straightforward method to differentiate between macromolecular crystals and salt crystals [1].

Protein and salt crystals differ substantially in their solvent content. Small crystal dyes, like JBS Deep Purple, are able to permeate the solvent channels of a protein, thus coloring the protein purple. In contrast, salt crystals are tightly packed and do not possess large solvent channels. They will therefore remain colourless.

Usage:

Simply add 0.5-1 µl of JBS Deep Purple to the crystallization drop containing the crystals of interest.

Coloring Time:

JBS Deep Purple colors protein crystals after a few minutes. Even if the color of the solution is only faintly purple under the microscope, proteins will be stained within 5-15 min.

Very occasionally, it has been reported that protein crystals did not absorb crystal dyes [2].

Selected References:

- [1] Wilkosz *et al.* (1995) Preliminary characterization of EcoRI-DNA co-crystals: incomplete factorial design of oligonucleotide sequences. *Acta Cryst. D* **51**:938.
 [2] Eckert *et al.* (2003) Crystallization and preliminary X-ray analysis of



JBS Deep Purple

Crystal Violet

Alicyclobacillus acidocaldarius endoglucanase CelA. *Acta Cryst. D* **59**:139.