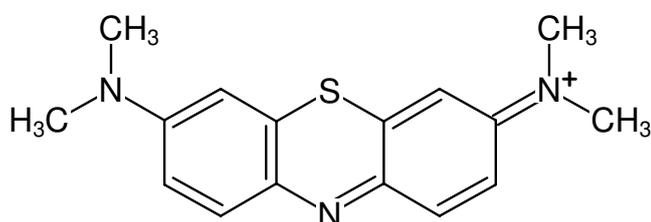


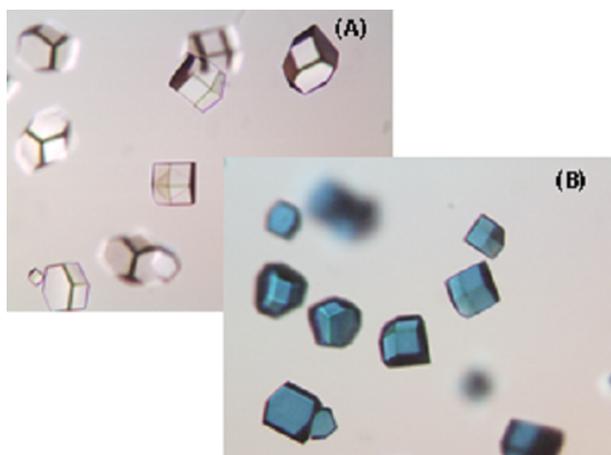
**JBS True Blue**

Methylene Blue hydrated

Cat. No.	Amount
CO-301	300 µl

Cl<sup>-</sup>

Structural formula of JBS True Blue



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:** C<sub>16</sub>H<sub>18</sub>ClN<sub>3</sub>S \* H<sub>2</sub>O**Molecular Weight:** 319.85 g/mol (anhydrous)**CAS#:** 61-73-4**EC number:** 200-515-2**Applications:**

JBS True Blue 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 True Blue, are able to permeate the solvent channels of a protein, thus coloring the protein blue. In contrast, salt crystals are tightly packed and do not possess large solvent channels. They will therefore remain colourless.

**Usage:**

Simply add 0.5 µl of JBS True Blue to the crystallization drop containing the crystals of interest.

**Coloring Time:**

JBS True Blue colors protein crystals after a few minutes. Even if the color of the solution is only faintly blue 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].

**Safety Information:**

Methylene Blue is a harmful substance if swallowed. Work carefully and wear gloves when handling the dye. Under fire conditions it emits toxic fumes.



## JBS True Blue

Methylene Blue hydrated

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