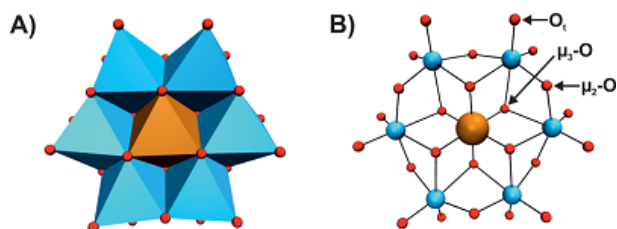




Anderson-Evans polyoxotungstate

Cat. No.	Amount
X-TEW-5	5 mg



Polyhedral (A) and ball and stick (B) representation of $[\text{TeW}_6\text{O}_{24}]^{6-}$ with different coordination modes of the oxygen atoms in panel B.^[1] Figure used by courtesy of Prof. Annette Rompel, University of Vienna.

For in vitro use only!

Shipping: shipped at ambient temperature

Storage Conditions: store at ambient temperature

Shelf Life: 12 months

Molecular Formula: $\text{Na}_6[\text{TeW}_6\text{O}_{24}] \times 22 \text{H}_2\text{O}$

Molecular Weight: 2148.90 g/mol

Solubility: 100 mM in water.

Dissolve 5 mg $\text{Na}_6[\text{TeW}_6\text{O}_{24}] \times 22 \text{H}_2\text{O}$ in 23 μl water to achieve a 100 mM stock solution.

Description:

The Anderson-Evans polyoxotungstate $[\text{TeW}_6\text{O}_{24}]^{6-}$ (TEW) is a universal and flexible additive for protein crystallization. With its planar structure and high negative charge, it has proven to promote crystal contacts, improve crystal quality and further provides a valuable anomalous signal for phasing due to 6 tungsten atoms^[1].

Related Products:

XP Up Screen, #CS-351

XP Screen, #CS-350

Selected References:

- [1] Bijelic *et al.* (2017) Ten Good Reasons for the Use of the Tellurium-Centered Anderson-Evans Polyoxotungstate in Protein Crystallography. *Acc. Chem. Res.* **50**:1441.
- [2] Molitor *et al.* (2016) Aurone synthase is a catechol oxidase with hydroxylase activity and provides insights into the mechanism of plant polyphenol oxidases. *Proc. Natl. Acad. Sci.* **113**:E1806.
- [3] Molitor *et al.* (2016) *In situ* formation of the first proteinogenically functionalized $[\text{TeW}_6\text{O}_{24}\text{O}_2(\text{Glu})]^{7-}$ structure reveals unprecedented chemical and geometrical features of the Anderson-type cluster. *Chem. Commun.* **52**:12286.
- [4] Molitor *et al.* (2015) Crystallization and preliminary crystallographic analysis of latent, active and recombinantly expressed aurone synthase, a polyphenol oxidase, from *Coreopsis grandiflora*. *Acta Cryst. F* **71**:746.
- [5] Bijelic *et al.* (2015) Hen Egg-White Lysozyme Crystallisation: Protein Stacking and Structure Stability Enhanced by a Tellurium(VI)-Centred Polyoxotungstate. *ChemBioChem* **16**:233.
- [6] Mauracher *et al.* (2014) Latent and active *ab*PP04 mushroom tyrosinase cocrystallized with hexatungstotellurate(VI) in a single crystal. *Acta Cryst. D* **70**:2301.
- [7] Mauracher *et al.* (2014) Crystallization and preliminary X-ray crystallographic analysis of latent isoform PPO4 mushroom (*Agaricus bisporus*) tyrosinase. *Acta Cryst. F* **70**:263.