

XP Screen

Crystallization Screen for improved Crystal Quality and Phasing

The XP Screen provides 96 of the most prominent crystallization conditions upgraded with the Anderson–Evans polyoxotungstate $[\text{TeW}_6\text{O}_{24}]^{6-}$ (TEW) at 1 mM final concentration. This additive has proven to promote crystal contacts, improve crystal quality and further provides a valuable anomalous signal for phasing due to 6 tungsten atoms. Optimized for TEW stability, the XP Screen improves crystallization and crystal diffraction quality of challenging targets.

Ordering Information

[Order Online »](#)

JBS-CS-350	1Kit
------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com