

Anderson-Evans polyoxotungstate

Ordering Information

[Order Online »](#)

JBS-X-TEW-5	5mg
-------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

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

MW 2148.9

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.

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