

# ATRX (human), (recombinant)

ATRX is a subunit of the ATRX:DAXX complex involved in H3.3 deposition, heterochromatin maintenance and transcriptional regulation. It belongs to a family of SNF2-related nucleosome remodeling enzymes (DNA-dependent ATPases). These ATP-dependent motor proteins are major factors of chromatin structure and transcriptional control in vivo. SNF2-related factors have been implicated in a variety of cancers in humans.

## Ordering Information

[Order Online »](#)

ENZ-PRT352-0050	50µl
-----------------	------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Handling</b>	Thaw and freeze quickly, in room-temperature water bath and liquid nitrogen, respectively; keep on ice when not in storage; split in small aliquots upon first thawing; avoid excessive freeze-thaw cycles.
<b>Long Term Storage</b>	-80°C
<b>Shipping</b>	Dry Ice

## Product Details

<b>Alternative Name</b>	Transcriptional regulator ATRX, X-linked helicase II, XH2, X-linked nuclear protein, XNP, Znf-HX
<b>Formulation</b>	Liquid. In 20 mM Tris-HCl, pH 7.9, 150 mM NaCl, 15% glycerol, 2 mM MgCl <sub>2</sub> , 0.2 mM EDTA, 1 mM DTT, 10 mM beta-glycerophosphate, 0.4 mM PMSF, 1 mM benzamidine and 0.4 mg/ml recombinant human insulin.
<b>MW</b>	~267 kDa
<b>Purity Detail</b>	Purified by immunoaffinity chromatography.
<b>Source</b>	Produced in insect (Sf9) cells. Human ATRX (aa 3-2375) fused to 6xHis and FLAG-tag at the N-terminus.
<b>UniProt ID</b>	P46100-3



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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