

EP400 (human), (recombinant)

EP400 is a subunit of the NuA4 histone acetyltransferase complex involved in nucleosome remodeling, core histone post-translational modifications 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-PRT347-0050	50µl
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	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.
Long Term Storage	-80°C
Shipping	Dry Ice

Product Details

Alternative Name	E1A-binding protein p400, CAG repeat protein 32, CAGH32, hDOM, KIAA1490, KIAA1818, p400 kDa SWI2/SNF2-related protein, Trinucleotide repeat-containing gene 12 protein, TNRC12
Formulation	Liquid. In 20 mM Tris-HCl, pH 7.9, 150 mM NaCl, 15% glycerol, 2 mM MgCl ₂ , 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.
MW	~333 kDa
Purity Detail	Purified by immunoaffinity chromatography.
Source	Produced in insect (Sf9) cells. Human EP400 (aa 2-3040) fused to 6xHis and FLAG-tag at the N-terminus.
UniProt ID	Q96L91-4



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)