

AMH (human), (recombinant)

AMH (anti-Müllerian hormone) is a member of the TGF- β family of genes and has been shown to be present in both males and females. It is known to be expressed in Sertoli cells and is involved in male sexual differentiation. The hormone is present in the ovary of females and has been shown to be a marker of female fertility. AMH was found to inhibit primordial follicle assembly and decrease the initial primordial follicle pool size. Women with polycystic ovarian syndrome (PCOS) exhibit elevated levels of Anti-Müllerian hormone (AMH) which has been found to desensitize ovarian follicles. It has also been shown to inhibit growth of cervical cancer cells via the MIS type II receptor.

Citations: 1

[View Online »](#)

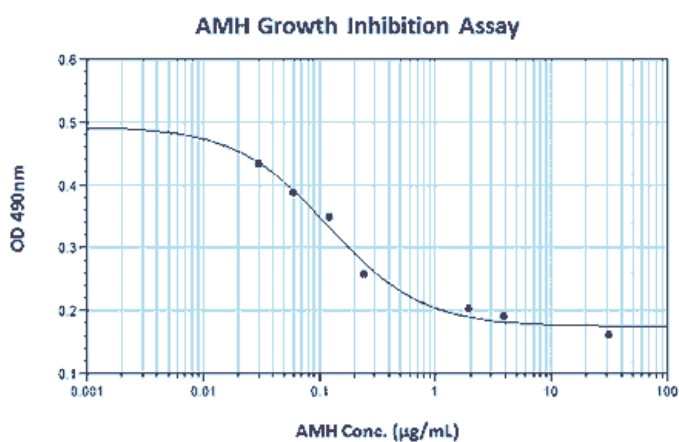
Ordering Information

[Order Online »](#)

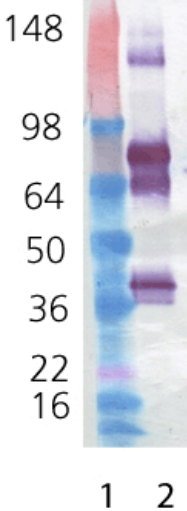
ENZ-PRT195-0010	10 μ g
-----------------	------------

Manuals, SDS & CofA

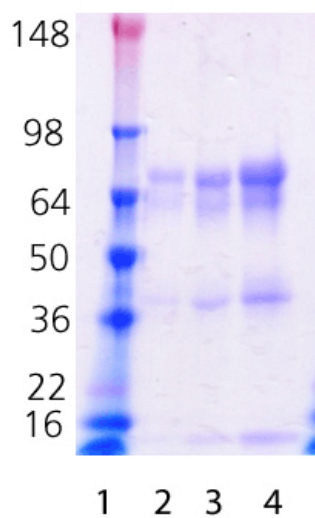
[View Online »](#)



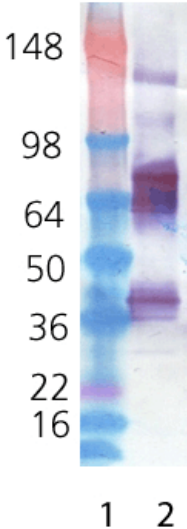
Caski cell growth inhibition assay after treatment with AMH (ENZ-PRT195): Caski cells were placed in stationary media at 5000 cells/well in 96 well plates. They were treated with a titrated concentration of AMH diluted in cell culture media. The cells were grown at 37°C / 5% CO₂ for 48 hours. Inhibition of growth was measured using a modified MTT cell viability assay. The ED₅₀ range is between 0.05 – 1.0 µg/mL.



Western blot analysis of AMH probed with mouse anti-HIS antibody. Lane 1: MWM, Lane 2: 100 ng AMH (Prod. No. ENZ-PRT195) probed HIS antibody.



Coomassie stained SDS-PAGE. Lane 1: Molecular weight marker, Lane 2: 1.0 µg, Lane 3: 2.0 µg, Lane 4: 5.0 µg AMH (Prod. No. ENZ-PRT195).



Western blot analysis of AMH probed with rabbit anti-FLAG antibody. Lane 1: MWM, Lane 2: 100 ng AMH (Prod. No. ENZ-PRT195) probed FLAG antibody.

Handling & Storage

Use/Stability	Stable for at least 6 months after receipt when stored at -20°C. After reconstitution, the protein is stable for 1 week at 4°C. For longer-term storage and/or applications sensitive to protein oligomerization, prepare aliquots and store at -80°C after reconstitution. Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	anti-Müllerian hormone, Mullerian-inhibiting substance
Biological Activity	Inhibits proliferation of Caski cells (ED ₅₀ range is between 0.05-1.0 µg/mL)
Concentration	0.1mg/ml when reconstituted as suggested.
Endotoxin Content	≤100EU/mg
Formulation	Lyophilized. Contains PBS.
MW	Expected is 61.2 kDa (main band is 68 kDa with minor bands at 60 kDa and 38 kDa)
Purity	≥90% (SDS-PAGE)
Purity Detail	Purified by multi-step chromatography.
Reconstitution	Reconstitute with 100µl sterile distilled water.
Source	Produced in HEK 293 cells. AMH (aa 19-560) is fused at the N-terminus to a FLAG [®] -tag and His-tag.
Technical Info / Product Notes	FLAG is a registered trademark of Sigma-Aldrich Co.
UniProt ID	P03971



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)