

ACTH monoclonal antibody (AH26)

This antibody specifically targets Synacthen (amino acids 1-24 of ACTH) and does not cross-react with CLIP (amino acids 17-39 of ACTH). It effectively stains corticotrophs in the adenohypophysis and proves valuable in the categorization of pituitary adenomas.

ACTH, or adrenocorticotrophic hormone, is a 39 amino acid active peptide hormone produced and secreted by the anterior pituitary gland. It plays a crucial role in the regulation of the adrenal glands and the synthesis of cortisol, a hormone involved in various physiological processes, including the body's response to stress. ACTH stimulates the adrenal cortex to release cortisol into the bloodstream. POMC (pro-opiomelanocortin or corticotropin-lipotropin) is a 267 amino acid polypeptide hormone precursor that goes through extensive, tissue-specific posttranslational processing by convertases. POMC is cleaved into ten hormone chains named NPP, ACTH, α -MSH (melanocyte stimulating hormone), β -MSH, γ -MSH, CLIP (corticotropin-like intermediary peptide), lipotropin- β , lipotropin- γ , β -endorphin and Met-enkephalin.

This antibody is covered by our [Worry-Free Guarantee](#).

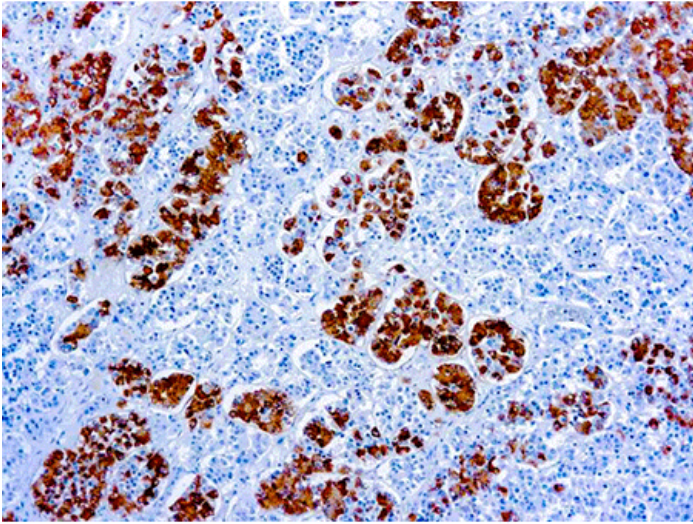
Ordering Information

[Order Online »](#)

ENZ-ABS851-1000	1ml
ENZ-ABS851-0100	100 μ l

Manuals, SDS & CofA

[View Online »](#)



Formalin fixed paraffin embedded human pituitary stained with Adrenocorticotrophic Hormone antibody.

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Adrenocorticotrophic hormone, Corticotropin
Application	IHC (PS)
Clone	AH26
Crossreactivity	Does not cross-react with CLIP (aa 17-39 of ACTH).
Formulation	This product is supplied as a purified immunoglobulin and contains less than 0.1% sodium azide as a preservative.
Host	Mouse
Immunogen	Synthetic peptide corresponding to amino acids 1-24 from the N-terminal of human ACTH.
Isotype	IgG1
Recommendation Dilutions/Conditions	For use with normal and neoplastic formalin-fixed, paraffin-embedded tissue sections. Recommended dilution 1:50-1:200.Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
Specificity	Specific to Synacthen (aa 1-24 of ACTH).
UniProt ID	P01189
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .



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)