



PRODUCT DATA SHEET

ACTH Recombinant monoclonal antibody (r57)

ENZ-ABS319

Product Number/Sizes

ENZ-ABS319-0100 100 µg

Product Details

ALTERNATIVE NAME: Adrenocorticotrophic hormone, Corticotropin
CLONE: r57
HOST: Mouse
ISOTYPE: IgG1
IMMUNOGEN: An N-terminal peptide (within aa 1-24) from the Synacthen/ACTH region of human POMC.
UNIPROT ID: P01189
SPECIES REACTIVITY: Human, Mouse, Rat
SPECIFICITY: This mAb is specific to Synacthen (aa 1-24 of ACTH); does not cross-react with CLIP (aa 17-39 of ACTH).
APPLICATIONS: IHC (PS)
RECOMMENDED DILUTIONS/CONDITIONS: Immunohistochemistry (formalin-fixed paraffin sections) (0.5-1 µg/mL) for 30 minutes at RT
Suggested dilutions/conditions may not be available for all applications.
Optimal conditions must be determined individually for each application.
PURITY DETAIL: Protein G affinity purified.
FORMULATION: Liquid. In 1X PBS containing 0.1 mg/ml BSA (US sourced) and 0.05% sodium azide.
SHIPPING: Blue Ice
LONG TERM STORAGE: +4°C
SCIENTIFIC BACKGROUND: ACTH (adrenocorticotrophic hormone, corticotropin) is a 39 amino acid active peptide produced by the anterior pituitary. 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. ACTH is also produced by cells of immune system (T-cells, B-cells, and macrophages) in response to stimuli associated with stress. Anti-ACTH is a useful marker in classification of pituitary tumors and the study of pituitary disease. It reacts with ACTH-producing cells (corticotrophs). It also may react with other tumors (e.g. some small cell carcinomas of the lung) causing paraneoplastic syndromes by secreting ACTH.
REGULATORY STATUS: RUO - Research Use Only

Revised 19-Dec-23

GLOBAL HEADQUARTERS

Enzo Life Sciences, Inc.
10 Executive Blvd
Farmingdale, NY 11735
USA
T 1-800-942-0430
T 1-631-694-7070
F 1-610-941-9252
E info-usa@enzolifesciences.com
www.enzolifesciences.com

EUROPE/ASIA

Enzo Life Sciences (ELS) AG
Industriestrasse 17, Postfach
CH-4415 Lausen
Switzerland
T +41/061 926 89 89
F +41/061 926 89 79
E info-ch@enzolifesciences.com
www.enzolifesciences.com