

Acidic Cytokeratin Recombinant monoclonal antibody (RMAK1-1)

ENZ-ABS335

Product Number/Sizes

ENZ-ABS335-0100 100 µg

Product Details

CLONE: RMAK1-1
HOST: Rabbit
ISOTYPE: IgGκ
IMMUNOGEN: An amino acid sequence common to acidic/type I keratins.
UNIPROT ID: Q7Z794
SPECIES REACTIVITY: Human
APPLICATIONS: IF, IHC (PS)
RECOMMENDED DILUTIONS/CONDITIONS: Immunofluorescence (0.5-1 µg/mL)
Immunohistochemistry (formalin-fixed paraffin sections) (0.5-1 µg/mL for 30 min at RT)
Suggested dilutions/conditions may not be available for all applications.
Optimal conditions must be determined individually for each application.
PURITY DETAIL: Protein A 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: There are two types of cytokeratins/keratins/CKs: the acidic type I cytokeratins and the basic or neutral type II cytokeratins. The subsets of cytokeratins which an epithelial cell expresses depends mainly on the type of epithelium, the moment in the course of terminal differentiation and the stage of development. Thus this specific keratin fingerprint allows the classification of all epithelia upon their keratin expression profile. Furthermore this applies also to the malignant counterparts of the epithelia (carcinomas), as the keratin profile tends to remain constant when an epithelium undergoes malignant transformation. The main clinical implication is that the study of the keratin profile by immunohistochemistry techniques is a tool of immense value widely used for tumor diagnosis and characterization in surgical pathology. [Wiki]
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