



# PRODUCT DATA SHEET

## Basic Cytokeratin Recombinant monoclonal antibody (KRTH/1576R)

ENZ-ABS331

### Product Number/Sizes

ENZ-ABS331-0100 100 µg

### Product Details

CLONE: KRTH/1576R  
HOST: Rabbit  
ISOTYPE: IgGκ  
IMMUNOGEN: Recombinant full length human KRT76.  
UNIPROT ID: Q01546  
SPECIES REACTIVITY: Human  
SPECIFICITY: This mAb recognizes basic (Type II or HMW) cytokeratins, which include 67kDa (CK1); 64kDa (CK3); 59kDa (CK4); 58kDa (CK5); 56kDa (CK6); 52kDa (CK8)  
APPLICATIONS: Flow Cytometry, IF, IHC (PS), WB  
RECOMMENDED DILUTIONS/CONDITIONS: Flow Cytometry (0.5-1 µg/10<sup>6</sup> cells in 0.1 mL)  
Immunofluorescence (1-2 µg/mL)  
Western blot (0.5-1 µg/mL)  
Immunohistochemistry (formalin-fixed paraffin sections) (0.25-0.5 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: Twenty human keratins are resolved with two-dimensional gel electrophoresis into acidic (pI 6.0) subfamilies. The acidic keratins have molecular weights (MW) of 56.5, 55, 51, 50, 50, 48, 46, 45, and 40kDa. Many studies have shown the usefulness of keratins as markers in cancer research and tumor diagnosis.  
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](mailto:info-usa@enzolifesciences.com)  
[www.enzolifesciences.com](http://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](mailto:info-ch@enzolifesciences.com)  
[www.enzolifesciences.com](http://www.enzolifesciences.com)