



PRODUCT DATA SHEET

Cadherin 16 Recombinant monoclonal antibody (CDH16/1532R)

ENZ-ABS360

Product Number/Sizes

ENZ-ABS360-0100 100 µg

Product Details

ALTERNATIVE NAME: Kidney-specific cadherin, Ksp-cadherin
CLONE: CDH16/1532R
HOST: Rabbit
ISOTYPE: IgGκ
IMMUNOGEN: Recombinant full length human Cadherin 16.
UNIPROT ID: O75309
SPECIES REACTIVITY: Human, Mouse, Rat
APPLICATIONS: Flow Cytometry, IF, IHC (PS)
RECOMMENDED DILUTIONS/CONDITIONS: Flow Cytometry (0.5-1 µg/10⁶ cells
Immunofluorescence (1-2 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: Cadherins form a superfamily of related glycoproteins that mediate calcium-dependent cell adhesion and transmit signals from the extracellular matrix to the cytoplasm. Cadherins have been implicated in embryogenesis, tissue morphogenesis, tissue structure maintenance, cell polarization, neoplastic invasiveness and metastasis, and membrane transport. It is suggested that CDH16 (kidney-specific cadherin, Cadherin 16) is a marker for terminal differentiation of the basolateral membranes of renal tubular epithelial cells. Within the kidney, CDH16 is found exclusively in the basolateral membrane of renal tubular epithelial cells and collecting duct cells, and not in glomeruli, renal interstitial cells, or blood vessels. CDH16 has been suggested to distinguish Chromophobe Renal-Cell Carcinoma from Oncocytoma.
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
T +41/061 926 89 79
E info-ch@enzolifesciences.com
www.enzolifesciences.com