p16^{INK4a} monoclonal antibody

Monoclonal antibody to p16 for use with IHC applications

p16INK4a (p16, CDKN2A) is a tumor suppressor that inhibits the progression of the cell cycle through the G1 phase. p16 expression is associated with high-risk human papillomavirus in cervical cancer, and head and neck tumors. It is frequently deleted or mutated in tumors such as melanomas, gliomas, esophageal, pancreatic, lung, urinary bladder carcinomas, and some types of leukemias; making p16 a key marker for cancer research.

This antibody is covered by our Worry-Free Guarantee.

Citations: 7

View Online »

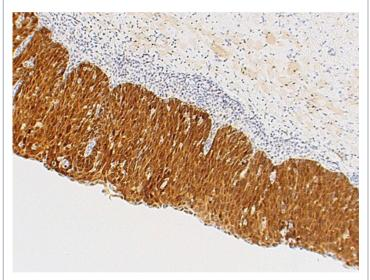
Ordering Information

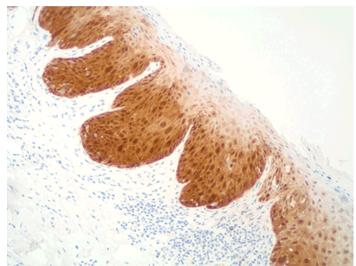
Order Online »

ENZ-ABS377-1000	1ml
ENZ-ABS377-0100	100μΙ

Manuals, SDS & CofA

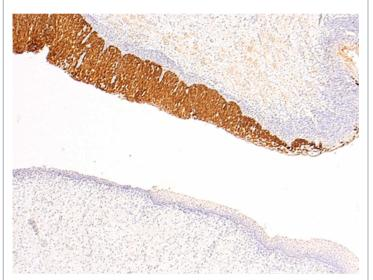
View Online »

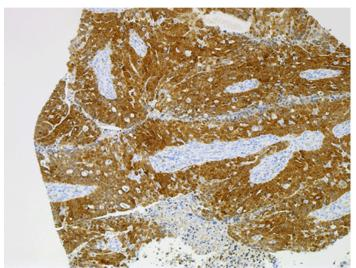




Formalin-fixed paraffin-embedded cervical tissue stained using Enzo's p16 antibody (ENZ-ABS377). IHC staining was performed on an automated Leica BOND III. Picture courtesy of Michael Priebs from Kantonsspital Aarau, Aarau, Switzerland.

Anal tissue incubated with p16INK4a monoclonal antibody and then stained with POLYVIEW® PLUS (anti-Mouse) HRP-DAB Kit (ENZ-KIT160).





Formalin-fixed paraffin-embedded cervical tissue stained using Enzo's p16 antibody (ENZ-ABS377). IHC staining was performed on an automated Leica BOND III. Picture courtesy of Michael Priebs from Kantonsspital Aarau, Aarau, Switzerland.

Formalin-fixed paraffin-embedded tissue punch from HPV-positive cervical carcinoma stained using Enzo's p16 antibody (ENZ-ABS377). IHC staining was performed on an automated Dako Autostainer Plus following antigen retrieval at high pH. Picture courtesy of Dr. Juana Maria Garcia Pedrero from Hospital Universitario Central de Asturias, Oviedo, Spain.

Handling & Storage

Handling Avoid freeze/thaw cycles.

Short Term Storage +4°C

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name p16, Cyclin-dependent kinase inhibitor 2A

Application IHC, WB

Application Notes For IHC applications: Dilute the concentrate as

appropriate in antibody diluent, and then apply 150 µL to

the slide, as indicated in the user's IHC protocol.

Formulation Liquid. In Tris buffer, pH 7.5, containing 1% BSA and

<0.1% sodium azide.

Host Mouse

Immunogen Recombinant p16.

lgG lgG

Purity Detail Affinity-purified.

Recommendation Dilutions/Conditions Immunohistochemistry (1:50-1:100)Optimal conditions

must be determined individually for each application.

Species Reactivity Human

UniProt ID P51480

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com