

# p16<sup>INK4a</sup> monoclonal antibody

Monoclonal antibody to p16 for use with IHC applications

p16<sup>INK4a</sup> (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.

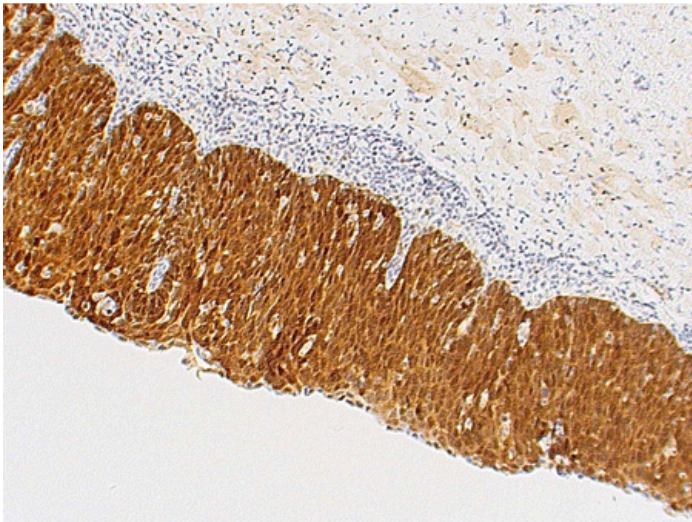
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 7 [View Online »](#)

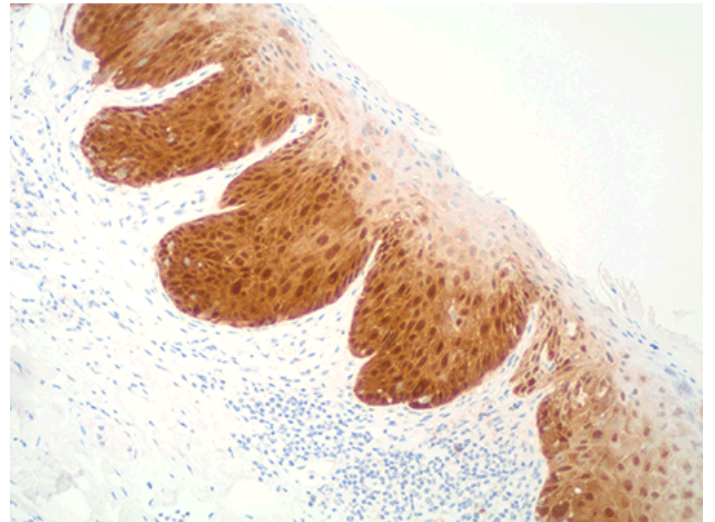
Ordering Information [Order Online »](#)

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

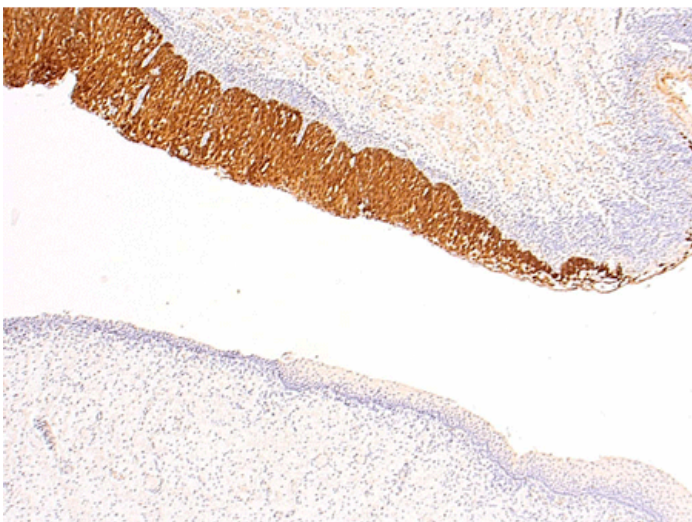
Manuals, SDS & CofA [View Online »](#)



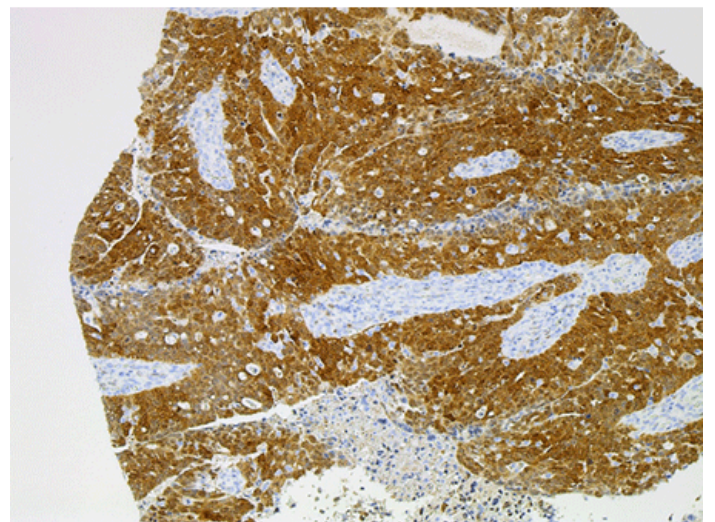
Formalin-fixed paraffin-embedded cervical tissue stained using Enzo's p16 antibody (ENZ-ABS377). IHC staining was performed on an automated Leica BOND III. Picture courtesy of Michael Priebs from Kantonsspital Aarau, Aarau, Switzerland.



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.

**Isotype** IgG

**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](#).



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)