

# p16<sup>INK4a</sup> monoclonal antibody (Ready-to-Use)

## Ready-to-use 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](#).

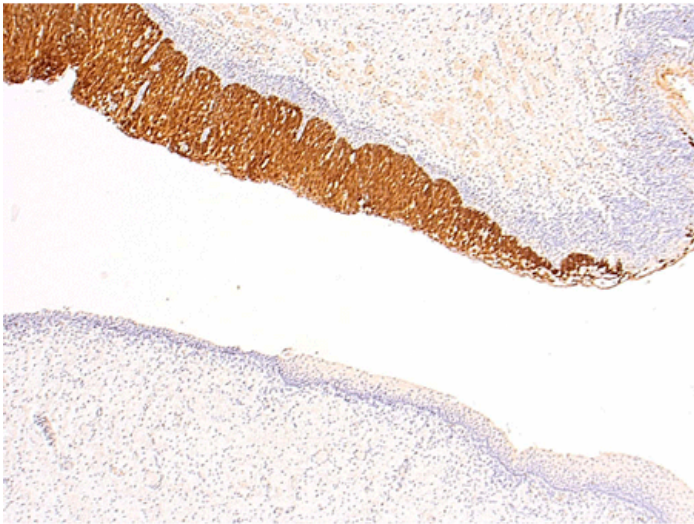
### Ordering Information

[Order Online »](#)

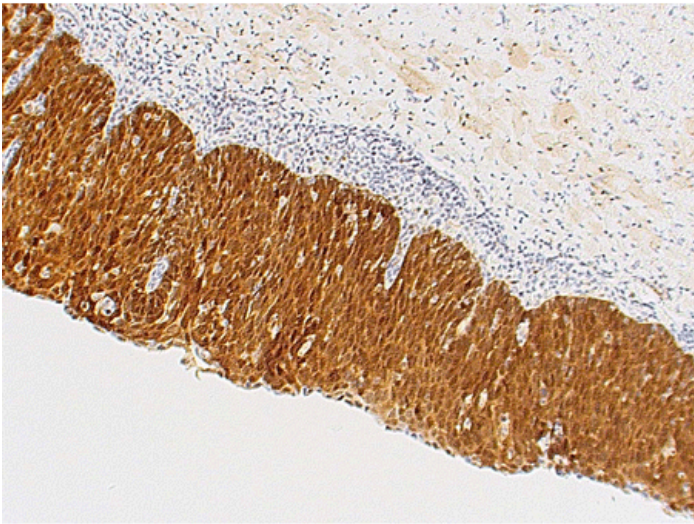
|                 |      |
|-----------------|------|
| ENZ-ABS664-0015 | 15ml |
|-----------------|------|

### Manuals, SDS & CofA

[View Online »](#)



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



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

## 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   |
| Formulation                         | Liquid. In Tris buffer, pH 7.5, containing 1% BSA and < 1% sodium azide.  |
| Host                                | Mouse   |
| Immunogen                           | Recombinant p16   |
| Isotype                             | IgG   |
| Purity Detail                       | Affinity purified.  |
| Recommendation Dilutions/Conditions | Immunohistochemistry (RTU)Suggested dilutions/conditions may not be available for all applications.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-  
fr@enzolifesciences.com](mailto:info-fr@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)

Last modified: May 29, 2024