PATHO-GENE[®] HPV type 16/18 probe

A mixture of biotin-labeled HPV 16 and HPV 18-specific probes for *in situ* hybridization

The PATHO-GENE® HPV type 16/18 probe is a mixture of biotin-labeled HPV 16 and HPV 18-specific probes in buffered formamide and hybridization enhancers.

Citations: 40

View Online »

Ordering Information

Order Online »

ENZ-32886	1ml
ENZ-32886-6000	6ml

Manuals, SDS & CofA

View Online »

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended. Stable for at least

one year after receipt when stored as recommended.

+4°C **Long Term Storage**

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Notes

The ENZO PATHO-GENE® HPV type 16/18 probe employs a mixture of the HPV 16 and HPV 18-specific probes that detect HPV types that have been associated with cervical intraepithelial neoplasia (CIN) and carcinoma in situ (CIS). The probe mixture is denatured and hybridized to the tissue section fixed and pretreated on the microscope slide. Specific hybridization between the HPV DNA probe and DNA in the specimen is determined by the detection of biotin.

Technical Info / Product Notes

HPV probe can be optimized for polymer detection system using diluent buffer, In situ hybridization buffer for HPV probes (Ready-to Use) (Prod. No. ENZ-33905).

This product or the use of this product is covered by one or more Enzo patents, including, but not limited to the following: U.S. Patent No. 4,994,373; Canadian Patent No. 1,309,672; and patents pending.

Last modified: May 29, 2024

