



PATHO-GENE® HPV Type 11 Probe

REF ENZ-GEN151-1000

INTENDED USE

Analyte Specific Reagent. Analytical and performance characteristics are not established.

REAGENTS PROVIDED

PATHO-GENE® HPV Type 11 probe is a biotin-labeled DNA probe designed to hybridize to HPV Type 11 DNA sequences. PATHO-GENE® HPV Type 11 Probe (0.1-2 µg/ml) is provided in buffer containing formamide, salts and dextran sulfate.

STORAGE & STABILITY

Store at 2°C to 8°C. When properly stored under these conditions, the product is stable to the expiration date indicated on the label. Do not use this product after the expiration date.

PRECAUTIONS



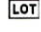




1. This product is an Analyte Specific Reagent (ASR). Analytical and performance characteristics are not established.
 2. The probe has affinity to HPV 11 genotype and may cross-react with known or uncharacterized viral types. It is the responsibility of the user to validate any test for its specific use.
 3. Specimens and all materials exposed to them should be considered potentially biohazardous, handled and disposed with proper precautions. Consult local and/or state authorities for the recommended method of disposal. Avoid contacting the skin and mucous membranes with reagents and specimens. Follow standard laboratory precautions to prevent exposure to eyes and skin. If reagents or specimens come in contact with sensitive areas, wash with copious amounts of water.
 4. Avoid microbial contamination of reagents as it may cause incorrect results.
 5. **Hazard-determining component of labeling:** Formamide
 - H360D May damage the unborn child.
 - P201 Obtain special instructions before use.
 - P202 Do not handle until all safety precautions have been read and understood.
 - P281 Use personal protective equipment as required.
 - P308+P313 IF exposed or concerned: Get medical advice/attention.
 - P405 Store locked up.
 - P501 This material and its container must be disposed of in a safe way.
- See also the product Safety Data Sheet (SDS) located at www.enzolifesciences.com.

INTELLECTUAL PROPERTY

PATHO-GENE® is a trademark of Enzo Biochem, Inc.

CONTACT INFORMATION

 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

Key to Symbols Used	
	Manufacturer
	Analyte Specific Reagent
	Batch code
	Catalogue number
	Use By
	Temperature Limitation
	Health Hazard