



## Human Papillomavirus 16<sup>GST</sup>

HPV 16  
recombinant, *E. coli*

Cat. No.	Amount
PR-1281	100 µg

### For *in vitro* use only!

**Shipping:** shipped on gel packs

**Storage Conditions:** store at -20 °C

**Additional Storage Conditions:** avoid freeze/thaw cycles

**Shelf Life:** 12 months

**Purity:** > 95 % (SDS-PAGE)

**Form:** liquid (Supplied in PBS, 100 mM K<sub>2</sub>CO<sub>3</sub> and 5 M urea)

### Description:

The Recombinant HPV-16 antigen is a full length protein expressed in *E. coli* having an MW of 56.7 kDa. The protein is fused to a 26 kDa GST-Tag.

Human papillomaviruses (HPVs) are a group family with more than 150 related viruses. HP 6 and 11 considered as low-risk HPVs can be sexually transmitted, causing warts to emerge on or around the genitals or anus, known as condylomata acuminata. HPV-6 major/large capsid antigen is used in clinical diagnosis to test the specific antibody to this virus induced by the virus infection, the positive reaction of this antibody is deemed as a marker for present and past infection. Furthermore, HPV-6 large capsid is used as a potential candidate for vaccine development.

### Amino Acid Sequence:

MQVTFIYILV ITCYENDVNV YHIFQMSLW LPSEATVYLP PVPVSKVVST  
 DEYVARTNIY YHAGTSRLLA VGHYPFIKK PNNNKILVPK VSGLQYRVFR  
 IHLDPNKFV FPDTSFYNDP TQRLVWACVG VEVGRGQPLG VGISGHPLLN  
 KLDDTENASA YAANAGVDNR ECISMDYKQT QLCLIGCKPP IGEHWGKGGSP  
 CTNVAVNP GD CPPELINTV IQDGDVHTG FGAMDFTTLQ ANKSEVPLDI  
 CTSICKYPDY IKMVSEPYGD SLFFYLRRREQ MFVRHLFNRA GTVGENVPDD  
 LYIKGSGSTA NLAASSNYFPT PSGSMVTS DA QIFNKPYWLQ RAQGHNNGIC  
 WGNQLFVTVV DTTRSTNMSL CAAISTSETT YKNTNFKEYL RHGEEYDLQFI  
 FQLCKITLTA DVMTYIHSMN STILEDWNFG LQPPPGTLE DTYRFVTQAI  
 ACQKHTPPAP KEDDPLKKYT FWEVNLKEKF SADLDQFPLG RKFLQAGLK  
 AKPKFTLGKR KATPTSSSTS TTAKRKRKL

### Selected References:

Munoz *et al.* (2010) Impact of human papillomavirus (HPV)-6/11/16/18 vaccine on all HPV-associated genital diseases in young women. *J Natl Cancer Inst.* **102**:325.