



Human Papillomavirus 16^{GST}

HPV 16
recombinant, *E. coli*

Cat. No.	Amount
PR-1281	100 µg

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Purity: > 90 % (SDS-PAGE)

Form: liquid (Supplied in PBS, 100 mM Arginine and 3 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 PVPVSKVST
 DEYVARTNIY YHAGTSRLLA VGHPYFPIKK PNNNKILVPK VSGLQYRVFR
 IHLDPNKFV FPDTSFYNDP TQRLVWACVG VEVGRGQPLG VGISGHPLLN
 KLDDTENASA YAANAGVDNR ECISMDYKQT QLCLIGCKPP IGEHWGKGGSP
 CTNVAVNPGD CPPELINTV IQDGDVMVHTG FGAMDFTTLQ ANKSEVPLDI
 CTSICKYPDY IKMVSEPYGD SLFFYLRRREQ MFVRHLFNRA GTVGENVPDD
 LYIKGSGSTA NCLASSNYFPT PSGSMVTSDA QIFNKPYWLQ RAQGHNNGIC
 WGNQLFVTWV DTTRSTNMSL CAAISTSETT YKNTNFKEYL RHGEEYDLQFI
 FQLCKITLTA DVMTYIHSMN STILEDWNFG LQPPPGGTLE DTYRFVTQAI
 ACQKHTPPAP KEDDPLKKYT FWEVNLKEKF SADLDQFPLG RKFLQAGLK
 AKPKFTLGKR KATPTTSSTS TTAKRKRKRL

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.