

**Human Papillomavirus 11<sup>GST</sup>**HPV 11  
recombinant, *E. coli*

Cat. No.	Amount
PR-1280	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:** > 95 % (SDS-PAGE)**Form:** liquid (Supplied in PBS, 3 M urea and 100 mM arginine)**Description:**

Recombinant HPV 11 antigen is a 58.1kDa protein covering the full length of HPV-11 major capsid, its N-terminus is fused with a GST tag. Human papillomaviruses (HPVs) are a group family with more than 150 related viruses. HPV 6 and 11 are considered as low-risk HPVs and 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 for 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.

**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.