

**Human Papillomavirus 6^{GST}**HPV 6
recombinant, *E. coli*

Cat. No.	Amount
PR-1279	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 50 mM arginine and 3M Urea)**Description:**

Recombinant HPV6 antigen is a 55.6 kDa protein covering the full length of HPV6 major capsid, its N-terminus is fused with a 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.