DATA SHEET





■ Human Papillomavirus 18 GST

HPV 18 recombinant, *E. coli*

Cat. No.	Amount
PR-1282	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 1x PBS and 2 mM Urea)

Description:

The recombinant HPV-18 (a 56kDa antigen) comprised from $\it E.~coli$ and fused to 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 acuminate. 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.

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.