

**HBVsAg adw2**

Recombinant Hepatitis B Surface Antigen adw subtype 2  
recombinant, *E. coli*

| Cat. No. | Amount |
|----------|--------|
| PR-1196  | 50 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:** > 90 % (SDS-PAGE)

**Form:** liquid (Supplied in 50 mM arginine and 1 M urea)

**Applications:**

Antigen in ELISA and Western blots, excellent antigen for detection of HBV with minimal specificity problems.

**Description:**

Recombinant HBVsAg adw antigen was mutated by replacing the Methionine residue at position 133 HBSAg is the surface antigen of the Hepatitis-B-Virus (HBV). The capsid of a virus has different surface proteins from the rest of the virus. The antigen is a protein that binds specifically on one of these surface proteins. It is commonly referred to as the Australian Antigen.

HbsAg subtype adw2 produced in *E. coli*, is a single, non-glycosylated, polypeptide chain containing 226 amino acids and having a molecular weight of approximately 23.0kDa. HBSAg adw2 migrates on SDS-PAGE as a 23kDa monomer band with minor amounts of dimer (46kDa) and trimer (69kDa) forms.

The HBSAg is fused to a 6 His tag on C-terminus and purified by proprietary chromatographic techniques.

**Specificity:** Immunoreactive with sera of HBV-infected individuals.