DATA SHEET





HBVsAg adw2

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

| Cat. No. | Amount |
|----------|--------|
| PR-1196 | 50 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 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 antigenof the Hepatitis-B-Virus (HBV). The capsidof 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.