

**HBsAg adw1 24 kDa**

Hepatitis B virus surface antigen adw1 subtype recombinant, *Pichia pastoris*

Cat. No.	Amount
PR-1402-S	200 µg
PR-1402-L	1 mg

For general laboratory use.

Shipping: shipped at 4 °C

Storage Conditions: store at 4 °C

Shelf Life: 6 months

Molecular Weight: about 24 kDa

Form: liquid (Supplied in Sodium Carbonate-Bicarbonate 20 mM pH 9.3, Sodium Chloride 100 mM, EDTA 3 mM, Glycerol 15 %)

Applications:

Suitable for use in ELISA. Each laboratory should determine an optimum working titer for use in its particular application. Other applications have not been tested but use in such assays should not necessarily be excluded.

Description:

The protein contains immunodominant region of HBV surface antigen adw1 subtype, amino acids 56-281.

Specificity:

Immunoreactive with sera of HBV-infected individuals