

**HBcore Ag 21 kDa**

Hepatitis B virus core antigen  
recombinant, *E. coli*

Cat. No.	Amount
PR-1400-S	200 µg
PR-1400-L	1 mg

**For in vitro use only!**

**Shipping:** shipped at 4 °C

**Storage Conditions:** store at 4 °C

**Shelf Life:** 6 months

**Molecular Weight:** about 21 kDa

**Form:** liquid (Supplied in Tris-HCl 50 mM pH 7.8, Sodium Chloride 50 mM, EDTA 0.5 mM, Glycerol 50 %)

**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 core antigen with 21 kDa, amino acids 2-149.

**Specificity:**

Immunoreactive with sera of HBV-infected individuals