

**HBeAg 44 kDa**

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

Cat. No.	Amount
PR-1401-S	200 µg
PR-1401-L	1 mg

For *in vitro* use only!

Shipping: shipped on blue ice

Storage Conditions: store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Molecular Weight: about 44 kDa

Purity: ≥ 85 %

Form: liquid (Supplied in Tris-HCl 50 mM pH 8.0, Sodium Chloride 60 mM, Glutathione 10 mM, Sarcosyl 0.25 %, 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 e antigen, amino acids 8-165. The protein has a GST tag located on the N terminus.

Specificity:

Immunoreactive with sera of HBV-infected individuals