

**TBEV gE 41 kDa**

Tick-borne encephalitis virus gE protein  
recombinant, *E. coli*

Cat. No.	Amount
PR-1450-S	200 µg
PR-1450-L	1 mg

**For general laboratory use.**

**Shipping:** shipped on gel packs

**Storage Conditions:** store at -20 °C

**Additional Storage Conditions:** avoid freeze/thaw cycles

**Shelf Life:** 6 months

**Molecular Weight:** about 41 kDa

**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 C-end immunodominant region of TBEV glycoprotein E, amino acids 296-414. The protein is fused to a GST tag.

**Specificity:**

Immunoreactive with sera of TBEV-infected individuals