

TBEV gE (residues 95-229)

Tick-borne Encephalitis Virus Glycoprotein E recombinant, *E. coli*

Cat. No.	Amount
PR-1227-1	1 mg

For *in vitro* use only
Quality guaranteed for 12 months
Store at -20°C

Avoid freeze / thaw cycles

Form

Liquid. Supplied in 20 mM MES pH 6.5, 8 M urea, 200 mM NaCl and 0.05% Tween-20.

Applications

Antigen in ELISA and Western blots, excellent antigen for detection of TBEV with minimal specificity problems.

Specificity

Immunoreactive with all sera of encephalitis virus-infected individuals.

Purity

>95% by SDS-PAGE

Description

The protein contains the Tick-borne Encephalitis Virus gE regions, amino acids: 95-229.

The protein is purified by proprietary chromatographic technique.

Background

Tick borne encephalitis virus (TBEV), is a human flavivirus causing tick-borne encephalitis (TBE), a viral infection of the central nervous system endemic in Europe and Asia. The recombinant E and C proteins triggered CD4⁺ but not CD8⁺ cells to proliferate and to produce IFN- γ and IL-5.

Selected References:

Gomez *et al.* (2003) Characterization of tick-borne encephalitis virusspecific human T lymphocyte responses by stimulation with structural TBEV proteins expressed in a recombinant baculovirus. *Viral Immunol.* **16**:407.