

**FSME (TBE) Virus Antigen (Premium)**

Strain Moscow B-4
native, Vero cell culture

Cat. No.	Amount
PR-BA112-S	250 µg
PR-BA112-L	1 mg

For *in vitro* use only!

Shipping: shipped on dry ice

Storage Conditions: store at -80 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Form: liquid

Culture system: Vero cell culture

Protein concentration: Lot specific. Determined by Bradford assay (Biorad) using BSA as standard.

Preservation: None

Analysis: ELISA

Preparation: Supernatant from infected culture is concentrated by centrifugation and subsequently purified by chromatography. The resulting virus preparation is presented in glycine buffer at pH 9.5. This preparation contains detergents.

Inactivation: This product has been treated with Beta-propiolactone. Confirmed inactive by attempted growth under original culture conditions. Since no procedure can guarantee absolute sterility, the reagent should be handled with appropriate safety precautions.