

**FSME (TBE) Virus Antigen**Strain Moscow B-4
native

| Cat. No. | Amount |
|-------------|--------|
| PR-BA112L-S | 250 µg |
| PR-BA112L-L | 1 mg |

For general laboratory use.**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. The resulting virus preparation is presented in glycine buffer at pH 9.5.**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.