

**Borrelia garinii** Antigen

native, Broth suspension

Cat. No.	Amount
PR-BA121GVS-S	250 µg
PR-BA121GVS-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:** Broth suspension**Protein concentration:** Lot specific. Determined by Bradford assay (Biorad) using BSA as standard.**Preservation:** None**Analysis:** ELISA**Preparation:** Harvested organisms are washed and detergent solubilised to produce a cytoplasmic extract enriched with lipoproteins and outer surface proteins (OSPs). The antigen is presented in phosphate buffered saline containing 1% n-octylglucoside.**Inactivation:** This product has been treated with detergent. 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.