

**Brucella abortus Antigen**S99 - Weybridge  
Fermentation

Cat. No.	Amount
PR-BA116-S	250 µg
PR-BA116-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 (Supplied in glycine buffer pH 9.5 containing n-octyl-β-D-glucopyranoside)**Culture system:** Fermentation**Protein concentration:** Lot specific. Determined by Bradford assay (Biorad) using BSA as standard.**Preservation:** None**Analysis:** ELISA**Preparation:** Bacterial lysate in glycine buffer (pH 9.5) containing n-octyl-β-D-glucopyranoside.**Inactivation:** This product has been treated with heat (1 hour, 65°C), phenol (0.5 % (w/w)) and 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.