

**Aspergillus fumigatus Antigen**

native, Liquid culture

Cat. No.	Amount
PR-BA132-S	250 µg
PR-BA132-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:** Liquid culture**Protein concentration:** Lot specific. Determined by Bradford assay (Biorad) using BSA as standard.**Preservation:** None**Analysis:** ELISA**Preparation:** Disrupted somatic material is clarified by centrifugation in phosphate buffered saline.**Inactivation:** Viable organisms are removed by centrifugation. 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.