

**Candida albicans Antigen**

wild strain  
native, Solid medium

Cat. No.	Amount
PR-BA117-S	250 µg
PR-BA117-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 (Supplied in 0.1 M glycine buffer pH 9.5)

**Culture system:** Solid medium

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

**Preservation:** None

**Analysis:** ELISA

**Preparation:** Organisms are washed, disrupted and subjected to an extraction process. The antigen consists of cytoplasmic and cell wall components, presented in alkaline pH glycine buffer.

**Inactivation:** This product has been inactivated by disruption. 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.