

**Chlamydia trachomatis**

EB & RB
native, Vero cell culture

Cat. No.	Amount
PR-BA1372-S	250 µg
PR-BA1372-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

CFT and SDS-PAGE

Preparation

Chlamydia trachomatis elementary and reticular bodies, purified from infected host cells by density gradient centrifugation. Presented in phosphate buffered saline pH 7.5.

Inactivation

Heat treatment and sonication. Inactivity is confirmed by attempted growth under original culture conditions. Since no procedure can guarantee absolute sterility, the reagent should be handled with appropriate safety precautions.

Agent

Chlamydia trachomatis LGV-II (434 strain)