

**Chlamydia MOMP 42 kDa**

Chlamydia trachomatis major outer membrane protein (MOMP) protein recombinant, *E. coli*

Cat. No.	Amount
PR-1451-S	200 µg
PR-1451-L	1 mg

For in vitro use only!

Shipping: shipped at 4 °C

Storage Conditions: store at 4 °C

Shelf Life: 6 months

Molecular Weight: about 42 kDa

Form: liquid (Supplied in Sodium Phosphate 100 mM pH 10.0, Tris-HCl 10 mM, Urea 8M, Imidasole 300mM)

Applications:

Suitable for use in ELISA. Each laboratory should determine an optimum working titer for use in its particular application. Other applications have not been tested but use in such assays should not necessarily be excluded.

Description:

The protein contains immunodominant region of Chlamydia trachomatis MOMP, amino acids 191-286. The protein is fused to a GST, 6 His, S-tag.

Specificity:

Immunoreactive with sera of Chlamydia trachomatis-infected individuals