

TRP p15

Treponema Pallidum Major Membrane Immunogen

Treponema pallidum, recombinant, *E. coli*

Cat. No.	Amount
PR-1239	100 µg

For *in vitro* use only
Quality guaranteed for 12 months
Store at -20°C

Avoid freeze / thaw cycles

Form

Liquid. Supplied in 10 mM Tris-HCl pH 8.0, 1 mM EDTA, 1 mM DTT and 8 M urea.

Application

Antigen in ELISA and Western blots, excellent antigen for detection of *T. pallidum* with minimal specificity problems.

Specificity

Immunoreactive with sera of *T. pallidum* infected individuals.

Purity

>90% by SDS-PAGE

Description

The protein contains the *Treponema pallidum* p15 immunodominant regions.

The protein is purified by proprietary chromatographic technique.

Background

Syphilis is a chronic, complex sexually transmitted disease of humans caused by the spirochetal bacterium *Treponema pallidum*. Humans are the only known reservoir for *T. pallidum*.

The 15 kDa lipoprotein of *Treponema pallidum* is a major immunogen during natural syphilis infection in humans and experimental infection in other hosts.

The humoral and cellular immune responses to this molecule appear late in infection as resistance to reinfection is developing.

Selected References:

- Centurion-Lara *et al.* (1997) Conservation of the 15-kilodalton lipoprotein among *Treponema pallidum* subspecies and strains and other pathogenic treponemes: genetic and antigenic analyses. *Infect. Immun.* **65**:1440.
- Baughn *et al.* (1996) Epitope mapping of B-cell determinants on the 15-kilodalton lipoprotein of *Treponema pallidum* (Tpp15) with synthetic peptides. *Infect. Immun.* **64**:2457.
- Fitzgerald (1992) Effects of cefetamet (Ro 15-8074) on *Treponema pallidum* and experimental syphilis. *Antimicrob. Agents Chemother.* **36**:598.
- Purcell *et al.* (1990) Lipid modification of the 15 kiloDalton major membrane immunogen of *Treponema pallidum*. *Mol. Microbiol.* **4**:1371.
- Purcell *et al.* (1989) Molecular cloning and characterization of the 15-kilodalton major immunogen of *Treponema pallidum*. *Infect. Immun.* **57**:3708.