

**TRP p15 partial^{His}**

Treponema Pallidum Major Membrane Immunogen
Treponema pallidum, recombinant, *E. coli*

Cat. No.	Amount
PR-1271	100 µg

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Molecular Weight: 48 kDa

Purity: > 90 % (SDS-PAGE)

Form: liquid (Supplied in 70 mM Tris-HCl pH 8.0, 50 mM NaCl, 50% glycerol and 1.5M urea)

Applications:

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

Description:

The *E. coli* derived recombinant 6xHis-tag fusion protein is a multimer having a molecular mass of 48kDa and containing the *T. pallidum* p15 immunodominant regions and six histidines fused at the C- terminus.

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.

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

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.