

**TRP p41 mosaic^{GST}**

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

Cat. No.	Amount
PR-1270	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

Purity: > 90 % (SDS-PAGE)

Form: liquid (Supplied in 25 mM Tris-HCl pH 8.0, 60 mM NaCl and 50% glycerol)

Applications:

Treponema Pallidum protein is suitable for ELISA and Western blots, excellent antigen for detection of *T. Pallidum* with minimal specificity problems.

Description:

The *E. coli* derived recombinant protein contains the outer membrane *T. Pallidum* p41 immunodominant regions. The protein is fused with a GST-tag.

Background: *Treponema pallidum* is a gram-negative spirochaete bacterium and is considered to be metabolically crippled. There are at least four known subspecies: *T. pallidum pallidum*, *T. pallidum pertenue*, *T. pallidum carateum* and *T. pallidum endemicum*. The helical structure of *T. pallidum pallidum* allows it to move in a corkscrew motion through viscous mediums such as mucus. *Treponema pallidum* sub sp. *pallidum* has one of the smallest bacterial genomes at 1.14 million base pairs (Mb) and has limited metabolic capabilities, reflecting its adaptation through genome reduction to the rich environment of mammalian tissue.

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

Selected References:

Poltavchenko *et al.* (2004) Dynamics of humoral immune response to *Treponema pallidum* proteins p17 and p41 at early stages of syphilis. *Zh. Mikrobiol. Epidemiol. Immunobiol.* 3:52.