

**Tp p47 47 kDa**

Treponema pallidum p47 protein
recombinant, *E. coli*

Cat. No.	Amount
PR-1428-S	200 µg
PR-1428-L	1 mg

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 6 months

Molecular Weight: about 47 kDa

Purity: 60 - 85 %

Form: liquid (Supplied in Sodium Phosphate 25 mM pH 8.0, Imidazole 150 mM, Sodium Chloride 150 mM, Glycerol 50 %)

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 *Treponema pallidum* p47, amino acids 33-375. The protein is fused to a 6 His TAG.

Specificity:

Immunoreactive with sera of *Treponema pallidum*-infected individuals