# **DATA SHEET**





# ■ Tp p47 164 kDa

Treponema pallidum p47 protein recombinant, *E. coli* 

| Cat. No.  | Amount |
|-----------|--------|
| PR-1423-S | 200 μg |
| PR-1423-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 164 kDa

**Purity:** 60 - 85 %

Form: liquid (Supplied in Tris-HCl 5 mM pH 8.0, Urea 4M, EDTA 0.5

mM, DTT 10 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  $\beta$ -galactosidase + 6 His tag.

# Specificity: