



P.O. Box 2746
Pinetop, AZ 85935
(Phone) 928-367-1200
(Fax) 928-367-1205

Catalog # 402P Antisera Pt GFP

Anti-*Ptilosarcus* GFP. Antibody to *Ptilosarcus* Green Fluorescent Protein

Background:

Ptilosarcus GFP is the 26 KDa green fluorescent protein cloned from the Sea Pen species known as *Ptilosarcus*. The purified recombinant protein was used as an antigen in conjunction with adjuvant to immunize Rabbits to high titer over several months.

Specificity:

By ELISA the sera can detect the *Ptilosarcus* GFP at a titer of 1:10,000.

Immunogen:

Purified *Ptilosarcus* GFP produced by recombinant means in *E. coli*.

Host:

Rabbit

Isotype:

IgG and IgM, mainly IgG species.

Species Reactivity:

Unknown cross reactivity.

Uses and Dilutions:

Western blotting: Use at a dilution of 1/1000. For best results, incubate membrane with diluted antibody in 5% w/v nonfat dry milk, 1X TBS, 0.1% Tween-20 at 4°C overnight with gentle agitation.

IHC(P): Use at a dilution of 1/2000. Optimal dilutions should be determined by the end user.

Positive Control:

Recombinant *Ptilosarcus* GFP.

Form:

250 microlitres

Concentration:

Unknown

Storage Buffer:

Rabbit sera

Storage:

Aliquot and store at -20 degrees Celcius. Avoid repeated freeze/thaw cycles.