



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

**Catalog # 403P Antisera Rr GFP**

Anti-*Renilla reniformis* GFP. Antibody to *Renilla reniformis* Green Fluorescent Protein

**Background:**

*R. reniformis* GFP is the 26 KDa green fluorescent protein cloned from the Atlantic Sea Pansy species known as *Renilla reniformis*. 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 *R. reniformis* GFP at a titer of 1:10,000.

**Immunogen:**

Purified *R. reniformis* 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 *R. reniformis* GFP.

**Form:**

250 microlitres

**Concentration:**

Unknown

**Storage Buffer:**

Rabbit sera

**Storage:**

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