



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

Catalog #401M Antisera Gluc

Anti-Gaussia Luciferase. Antibody to *Gaussia princeps* Luciferase

Background:

Gaussia Luciferase is the 19.9 KDa luciferase cloned from the deep sea copepod, *Gaussia princeps*.

The recombinant protein was used as an antigen in conjunction with adjuvant to immunize mice.

Specificity:

ELISA the sera can detect the Gaussia Luciferase at a titer of 1:10,000.

Immunogen:

Purified Gaussia Luciferase produced by recombinant means in bacteria.

Host:

mouse

Isotype:

IgG Kappa

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 2% w/v nonfat dry milk, 1X TBS, 0.1% Tween-20 at Room temperature for 1 hr.

IHC(P): Use at a dilution of 1/200 to 1:400, Optimal dilutions should be determined by the end user.

Positive Control:

Recombinant Gaussia Luciferase Proteins.

Form:

250 microlitres

Concentration:

Unknown

Storage Buffer:

DMEM

Storage:

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