

# PAb Control

Normal Rat IgG Control

Catalog code: pab-sctr

<https://www.invivogen.com/pab-control>

For research use only, not for diagnostic or therapeutic use

Version 21D20-MM

## PRODUCT INFORMATION

### Contents

- 200 µg purified rat IgG, provided azide-free and lyophilized.

### Formulation

H<sub>2</sub>O with 250 U/ml Pen and 250 µg/ml Strep

### Antibody resuspension

Add 1 ml of sterile phosphate-buffered saline (PBS) to obtain a concentration of 0.2 mg/ml.

### Storage and stability

- Product is shipped at room temperature. Lyophilized product should be stored at -20°C. Lyophilized product is stable for 1 year when properly stored.
- Resuspended product should be stored at -20°C for 1 year.

### Quality control

- Absence of neutralization activity has been tested using cellular assays.
- The absence of bacterial contamination (e.g. lipoproteins and endotoxins) has been confirmed using HEK-Blue™ TLR2 and HEK-Blue™ TLR4 cells.

## DESCRIPTION

Sera from Wistar rats were harvested and the IgG fraction purified by Protein G affinity chromatography. PAb Control is the control for use with PAb-hTLR2, PAb-hTLR4, and PAb-hTLR6. PAb Control was obtained from rats housed under the same conditions as the immunized rats used to produce the polyclonal TLR antibodies.

## APPLICATION

PAb Control is provided as a negative control for use in conjunction with InvivoGen's rat polyclonal antibodies. This control product should be used at the same concentration as the specific polyclonal antibody. Optimal dilutions should be determined for each application.

## RELATED PRODUCTS

Product	Description	Catalog Code
PAb-hTLR2	Human TLR2 Neutralizing antibody, Polyclonal Rat IgG	pab-htlr2
PAb-hTLR4	Human TLR4 Neutralizing antibody, Polyclonal Rat IgG	pab-htlr4
PAb-hTLR6	Human TLR6 Neutralizing antibody, polyclonal rat IgG	pab-htlr6

### TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Asia: +852 3-622-34-80

E-mail: [info@invivogen.com](mailto:info@invivogen.com)

