

PAG2 Polyclonal Antibody

(Catalog # A69915)

Background

PAG2 or a processed derivative of this molecule might represent a factor that binds the LH receptor.

Description

PAG2 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, pH 7.4.

Specificity

Bovine

Isotype

IgG

Uniprot ID

Q28057

Purification

>95%, Protein G purified

Immunogen

Recombinant Bovine Pregnancy-associated glycoprotein 2 protein (22-376AA)

Storage

Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Pregnancy-associated glycoprotein 2, PAG 2, PAG2

Application

ELISA
