
Immunoglobulin G2a Polyclonal Antibody

(Catalog # A50221)

Background

Immunoglobulin G (IgG) is a glycoprotein antibody that regulates immune responses such as phagocytosis and is also involved in the development of autoimmune diseases. Mouse IgGs have four distinct isotypes, namely, IgG1, IgG2a, IgG2b, and IgG3. IgG1 regulates complement fixation in mice. Mouse IgG2a antibodies react with anti-mouse whole serum and anti-mouse IgG2a.

Description

Immunoglobulin G2a Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Mouse

Isotype

IgG

Purification

Protein G purified

Immunogen

IgG2a purified from Mouse serum (Native Protein)

Storage

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

Alternative Names

Immunoglobulin G2a

Application

ELISA; recommended dilution: ELISA 1:1000-1:5000
