

rgpA Polyclonal Antibody

(Catalog # A56206)

Background

Thiol protease which is believed to participate in intracellular degradation and turnover of proteins. Its proteolytic activity is a major factor in both periodontal tissue destruction and in bacterial host defense mechanisms. Activates complement C3 and C5.

Description

rgpA Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Porphyromonas gingivalis

Isotype

IgG

Uniprot ID

P28784

Purification

>95%, Protein G purified

Immunogen

Recombinant Porphyromonas gingivalis Gingipain R1 protein (228-720AA)

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

Arg-gingipain, Gingipain 1, RGP-1, rgp1, rgpA

Application

ELISA