
GLA Polyclonal Antibody, HRP Conjugated

(Catalog #A53219)

Background

Alpha-galactosidase is a glycoside hydrolase enzyme that hydrolyses the terminal alpha-galactosyl moieties from glycolipids and glycoproteins.

Description

GLA Polyclonal Antibody, HRP Conjugated. HRP. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

P06280

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Alpha-galactosidase A protein (32-429AA)

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

Alpha-D-galactosidase A, Alpha-D-galactoside galactohydrolase, Melibiase, INN=Agalsidase, GLA

Application

ELISA; Recommended dilution: ELISA 1:500 - 1:1000
