
Vitamin D-Binding Protein Polyclonal Antibody, Biotin Conjugated

(Catalog # A57869)

Background

Multifunctional protein found in plasma, ascitic fluid, cerebrospinal fluid, and urine and on the surface of many cell types. In plasma, it carries the vitamin D sterols and prevents polymerization of actin by binding its monomers. DBP associates with membrane-bound immunoglobulin on the surface of B-lymphocytes and with IgG Fc receptor on the membranes of T-lymphocytes.

Description

Vitamin D-Binding Protein Polyclonal Antibody, Biotin Conjugated. Biotin. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

P02774

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Vitamin D-binding protein (19-474AA)

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

Gc-globulin, Group-specific component, GC

Application

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