
FABP3 Polyclonal Antibody, HRP Conjugated

(Catalog # A54251)

Description

FABP3 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

P05413

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Fatty acid-binding protein, heart protein (2-133AA)

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

Fatty acid-binding protein 3, Heart-type fatty acid-binding protein, H-FABP, Mammary-derived growth inhibitor, MDGI, Muscle fatty acid-binding protein, M-FABP, FABP3, FABP11, MDGI

Application

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