
TSPAN8 Polyclonal Antibody, Biotin Conjugated

(Catalog # A63727)

Background

extracellular exosome, integral component of plasma membrane, cell surface receptor signaling pathway

Description

TSPAN8 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

P19075

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Tetraspanin-8 protein (110-205AA)

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

Tetraspanin-8, Tspan-8, Transmembrane 4 superfamily member 3, Tumor-associated antigen CO-029, TSPAN8, TM4SF3

Application

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