
MUC16 Polyclonal Antibody, FITC Conjugated

(Catalog # A54127)

Description

MUC16 Polyclonal Antibody, FITC Conjugated. FITC. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

Q8WXI7

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Mucin-16 protein (21712-21973AA)

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

MUC-16, Ovarian cancer-related tumor marker CA125, MUC16, CA125

Application

ELISA; Recommended dilution: ELISA 1:100-1:500
