
TGFA Polyclonal Antibody, FITC Conjugated

(Catalog # A54055)

Description

TGFA 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

P01135

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Protransforming growth factor alpha protein (40-89AA)

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

EGF-like TGF, ETGF, TGF type 1

Application

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