
TGFB3 Polyclonal Antibody

(Catalog #A70903)

Background

This gene encodes a member of the TGF-beta family of proteins. The encoded protein is secreted and is involved in embryogenesis and cell differentiation. Defects in this gene are a cause of familial arrhythmogenic right ventricular dysplasia 1.

Description

TGFB3 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P10600

Purification

Affinity Purified

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 25-300 of human TGFB3 (NP_003230.1).

Storage

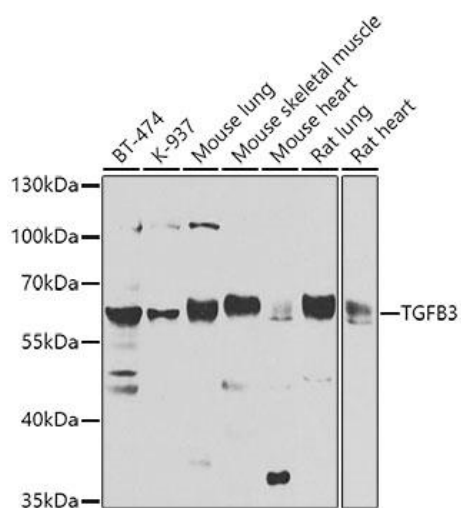
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

TGFB3; ARVD; ARVD1; LDS5; RNHF; TGF-beta3; transforming growth factor beta-3

Application

ELISA, WB, IHC: Recommended dilution; WB 1:500 - 1:2000, IHC 1:50-1:200



Western blot analysis of extracts of various cell lines, using TGFβ3 Polyclonal Antibody at 1:1000 dilution.

Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.

Lysates/proteins: 25ug per lane.

Blocking buffer: 3% nonfat dry milk in TBST.

Exposure time: 90s.