
RARG Polyclonal Antibody, FITC Conjugated

(Catalog A68677)

Background

Receptor for retinoic acid. Retinoic acid receptors bind as heterodimers to their target response elements in response to their ligands, all-trans or 9-cis retinoic acid, and regulate gene expression in various biological processes. The RAR/RXR heterodimers bind to the retinoic acid response elements (RARE) composed of tandem 5'-AGGTCA-3' sites known as DR1-DR5. In the absence of ligand, acts mainly as an activator of gene expression due to weak binding to corepressors. Required for limb bud development. In concert with RARA or RARB, required for skeletal growth, matrix homeostasis and growth plate function.

Description

RARG 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

P13631

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Retinoic acid receptor gamma protein (13-87AA)

Storage

Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Retinoic acid receptor gamma, RAR-gamma, Nuclear receptor subfamily 1 group B member 3, RARG, NR1B3

Application

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