
GADD45B Polyclonal Antibody

(Catalog #A62626)

Background

Involved in the regulation of growth and apoptosis. Mediates activation of stress-responsive MTK1/MEKK4 MAPKKK.

Description

GADD45B Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

O75293

Purification

Protein G purified

Immunogen

Recombinant Human Growth arrest and DNA damage-inducible protein GADD45 beta protein (1-160AA)

Storage

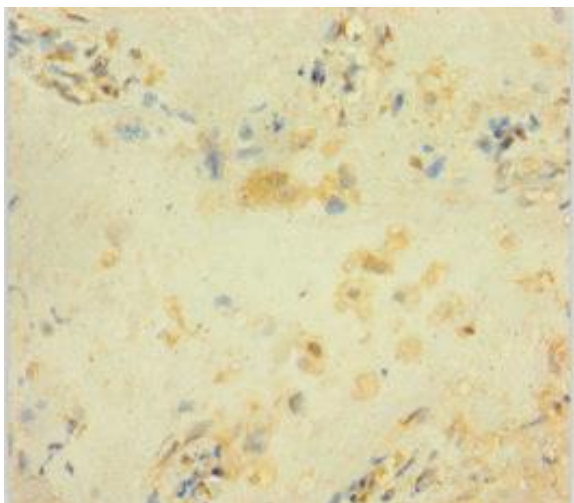
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

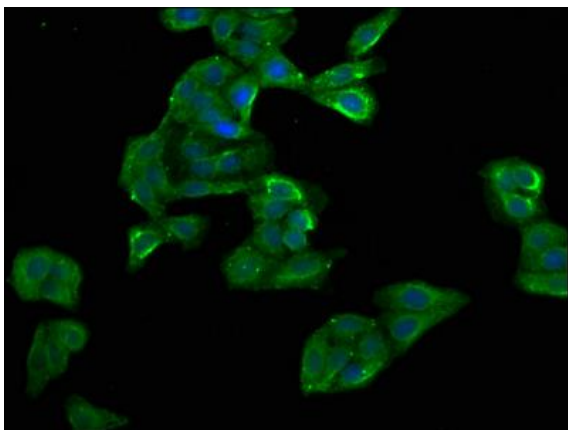
Growth arrest and DNA damage-inducible protein GADD45 beta, Myeloid differentiation primary response protein MyD118, Negative growth regulatory protein MyD118, GADD45B, MYD118

Application

ELISA, IHC, IF; Recommended dilution: IHC:1:20-1:200, IF:1:50-1:200



Immunohistochemistry of paraffin-embedded human placenta tissue using GADD45B Antibody at dilution of 1:100



Immunofluorescence staining of HepG2 cells with GADD45B Antibody at 1:100, counter-stained with DAPI. The cells were fixed in 4% formaldehyde, permeabilized using 0.2% Triton X-100 and blocked in 10% normal Goat Serum. The cells were then incubated with the antibody overnight at 4°C. The secondary antibody was Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L).