

HAS1 Polyclonal Antibody

(Catalog # A68551)

Background

Catalyzes the addition of GlcNAc or GlcUA monosaccharides to the nascent hyaluronan polymer. Therefore, it is essential to hyaluronan synthesis a major component of most extracellular matrices that has a structural role in tissues architectures and regulates cell adhesion, migration and differentiation. This is one of the isozymes catalyzing that reaction. Also able to catalyze the synthesis of chito-oligosaccharide depending on the substrate.

Description

HAS1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

Q92839

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Hyaluronan synthase 1 protein (151-271AA)

Storage

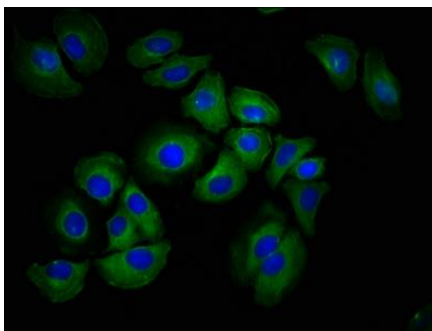
Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Hyaluronan synthase 1, Hyaluronate synthase 1, Hyaluronic acid synthase 1, HA synthase 1, HuHAS1, HAS1, HAS

Application

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



Immunofluorescence staining of A549 cells with HAS1 Polyclonal 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).