

AGT Polyclonal Antibody

(Catalog # A70590)

Background

The protein encoded by this gene, pre-angiotensinogen or angiotensinogen precursor, is expressed in the liver and is cleaved by the enzyme renin in response to lowered blood pressure. The resulting product, angiotensin I, is then cleaved by angiotensin converting enzyme (ACE) to generate the physiologically active enzyme angiotensin II. The protein is involved in maintaining blood pressure and in the pathogenesis of essential hypertension and preeclampsia. Mutations in this gene are associated with susceptibility to essential hypertension, and can cause renal tubular dysgenesis, a severe disorder of renal tubular development. Defects in this gene have also been associated with non-familial structural atrial fibrillation, and inflammatory bowel disease.

Description

AGT 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

P01019

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 35-285 of human AGT (NP_000020.1).

Storage

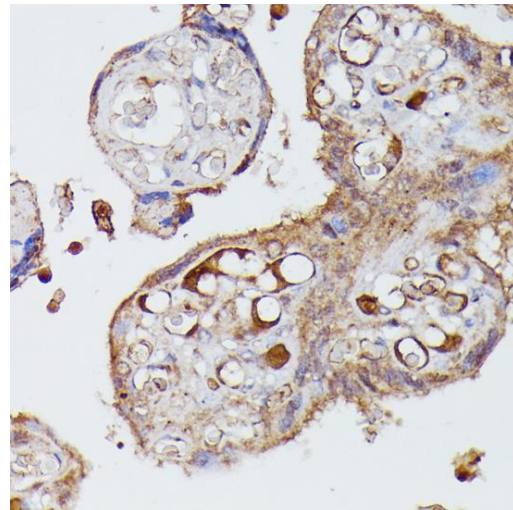
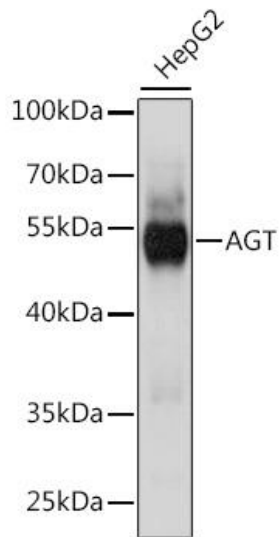
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

AGT; ANHU; SERPINA8; hFLT1; angiotensinogen

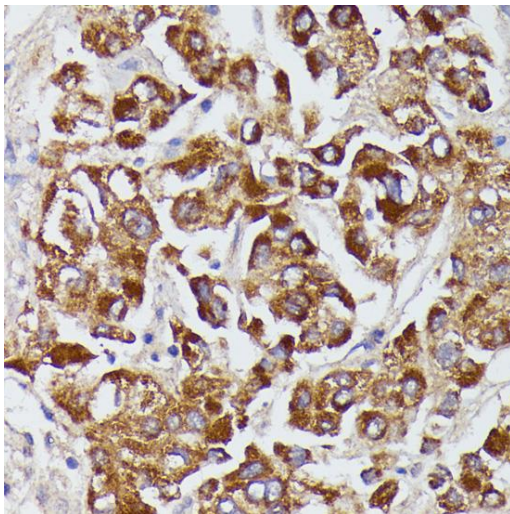
Application

WB, IHC, IF; Recommended dilution: WB 1:500 - 1:2000, IHC 1:50 - 1:100, IF 1:50 - 1:100



Immunohistochemistry of paraffin-embedded human placenta using AGT Polyclonal Antibody at dilution of 1:100 (40x lens).

Western blot analysis of extracts of HepG2 cells, using AGT 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: 30s.



Immunohistochemistry of paraffin-embedded human liver cancer using AGT Polyclonal Antibody at dilution of 1:100 (40x lens).