
Avpr1a Polyclonal Antibody

(Catalog # A53272)

Background

Receptor for arginine vasopressin. The activity of this receptor is mediated by G proteins which activate a phosphatidyl-inositol-calcium second messenger system. Involved in social memory formation.

Description

Avpr1a Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Preservative: 0.1mol/L NaCl, 15mmol/L NaN₃, PH7.2.

Specificity

Rat, Mouse

Isotype

IgG

Uniprot ID

P30560

Purification

Protein G purified

Immunogen

Recombinant Rat Vasopressin V1a receptor protein (7-52AA)

Storage

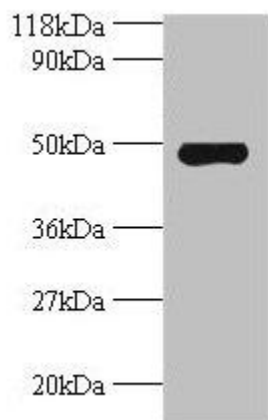
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

AVPR V1a Antidiuretic hormone receptor 1a Vascular/hepatic-type arginine vasopressin receptor Avpr1a

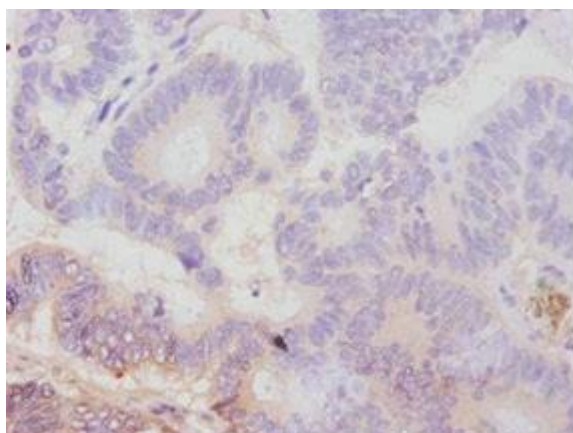
Application

ELISA, WB, IHC; Recommended dilution: WB:1:1000-1:5000, IHC:1:20-1:200



Western blot

All lanes: Avpr1a antibody at 2ug/ml + Mouse spleen tissue
Secondary
Goat polyclonal to rabbit IgG at 1/10000 dilution
Predicted band size: 48 kDa
Observed band size: 48 kDa



Immunohistochemistry of paraffin-embedded human colon cancer using Avpr1a Antibody at dilution of 1:100