
CHMP2A Polyclonal Antibody

(Catalog # A51805)

Background

Probable core component of the endosomal sorting required for transport complex III (ESCRT-III) which is involved in multivesicular bodies (MVBs) formation and sorting of endosomal cargo proteins into MVBs. MVBs contain intraluminal vesicles (ILVs) that are generated by invagination and scission from the limiting membrane of the endosome and mostly are delivered to lysosomes enabling degradation of membrane proteins, such as stimulated growth factor receptors, lysosomal enzymes and lipids. The MVB pathway appears to require the sequential function of ESCRT-O, -I, -II and -III complexes. ESCRT-III proteins mostly dissociate from the invaginating membrane before the ILV is released. The ESCRT machinery also functions in topologically equivalent membrane fission events, such as the terminal stages of cytokinesis and the budding of enveloped viruses (HIV-1 and other lentiviruses). ESCRT-III proteins are believed to mediate the necessary vesicle extrusion and/or membrane fission activities, possibly in conjunction with the AAA ATPase VPS4. Involved in HIV-1 p6- and p9-dependent virus release.

Description

CHMP2A 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

O43633

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Charged multivesicular body protein 2a protein (1-222AA)

Storage

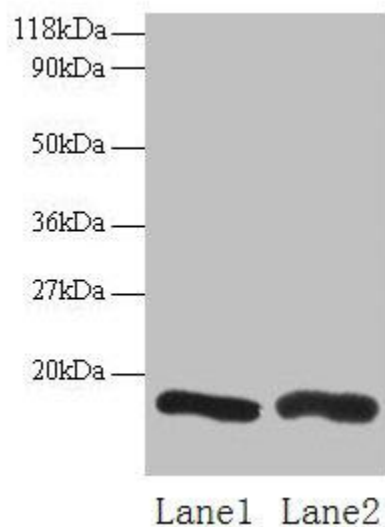
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

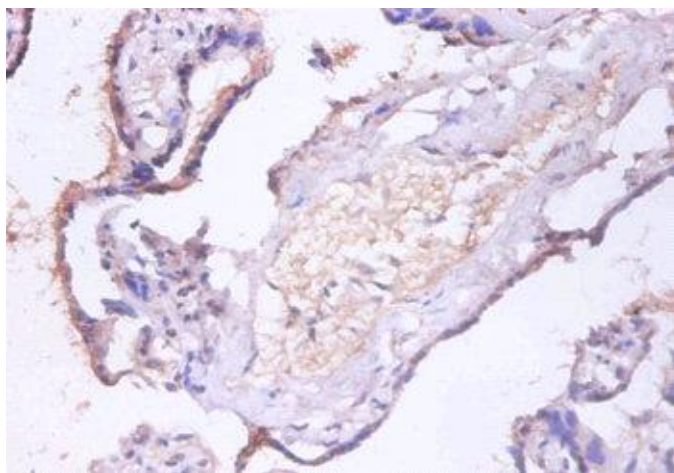
Charged multivesicular body protein 2a (Chromatin-modifying protein 2a) (CHMP2a) (Putative breast adenocarcinoma marker BC-2) (Vacuolar protein sorting-associated protein 2-1) (Vps2-1) (hVps2-1), CHMP2A, BC2 CHMP2

Application

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



All lanes: CHMP2A Polyclonal Antibody at 2ug/ml
Lane 1: EC109 whole cell lysate
Lane 2: 293T whole cell lysate
Secondary
Goat polyclonal to Rabbit IgG at 1/15000 dilution
Predicted band size: 24 kDa
Observed band size: 16kDa



Immunohistochemistry of paraffin-embedded human placenta at dilution of 1:100