
L1 Polyclonal Antibody

(Catalog # A52150)

Background

Forms an icosahedral capsid with a T=7 symmetry and a 50 nm diameter. The capsid is composed of 72 pentamers linked to each other by disulfide bonds and associated with L2 proteins. Binds to heparan sulfate proteoglycans on the basement membrane to provide initial virion attachment to target cells. Basement membrane is exposed only after epithelium trauma. Additionally, the alpha6 integrin complexed with either beta1 or beta4 integrin has been proposed to act as a coreceptor recognized by L1. Once attached, integrin complexed with beta4 integrin has been proposed to act as a coreceptor recognized by L1. Once attached, the virion enters the host cell via clathrin-mediated endocytosis and the genomic DNA is released to the host nucleus. The virion assembly takes place within the cell nucleus. Encapsulates the genomic DNA together with protein L2.

Description

L1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human papillomavirus type 18

Isotype

IgG

Uniprot ID

P06794

Purification

Protein G purified

Immunogen

Recombinant HPV18 Major capsid protein L1 protein (1-568AA)

Storage

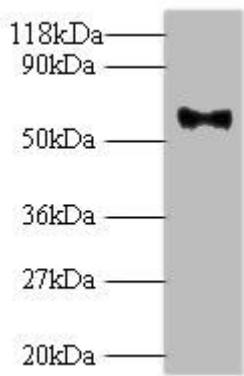
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

L1

Application

ELISA, WB; Recommended dilution: WB:1:1000-1:5000



Western blot

All lanes: L1 Polyclonal Antibody at 2ug/ml+293T whole cell lysate

Secondary

Goat polyclonal to Rabbit IgG at 1/10000 dilution

Predicted band size: 62kDa

Observed band size: 62kDa