

SLK Monoclonal Antibody [3E9]

[Catalog No. A-7168-100]

Description

Mouse monoclonal antibody raised against a partial recombinant SLK, clone 3E9.

Size

100 µg

Purification

Protein A purified

Immunogen

SLK (NP_055535, 1126 a.a. ~ 1236 a.a) partial recombinant protein with GST tag. MW of the GST tag alone is 26 KDa.

Cross Reactivity

Human

Isotype

IgG1 Kappa

Formulation

In 1x PBS, pH 7.2

Storage

Aliquot and store at -20°C or lower. Avoid repeated freezing and thawing. Multiple freeze/thaw cycles may result in decreased performance. Non-hazardous. No MSDS necessary.

Handling Recommendations

For maximum recovery of the products, centrifuge the vial prior to opening the cap.

Applications & Suggested Dilutions*

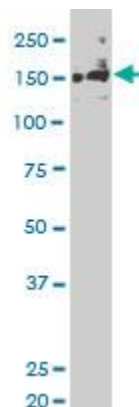
Western Blot: 0.5-1 µg/ml

ELISA: 0.25-1 µg/ml

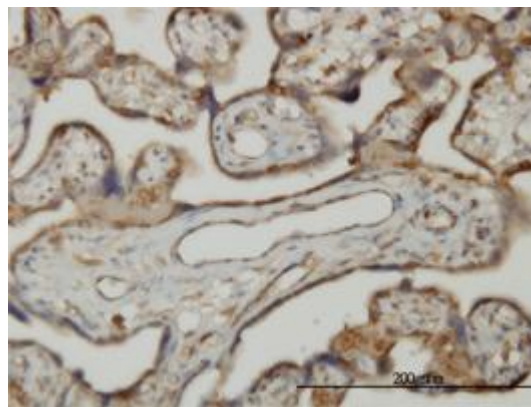
Immunohistochemistry (IHC): 2.5 µg/ml

**The end user is responsible for determining optimal working dilutions.*

Data



SLK Monoclonal Antibody, clone 3E9. Western Blot analysis of SLK expression in HeLa NE.



Immunoperoxidase of monoclonal antibody to SLK on formalin-fixed paraffin-embedded human placenta. [Antibody Concentration: 3 µg/ml]

This product is for research purposes only. Not intended for use in diagnostic procedures.