

Anti-NOTCH 4

Rabbit Polyclonal Antibody

Catalog # N44-363R

Lot # Z2014-20

Cited Applications

ELISA, IHC

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the NOTCH 4 protein

Cross Reactivity

Human

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

Rabbit polyclonal NOTCH4 antibody was raised against a peptide from the C terminal residues of human NOTCH4.

Formulation

PBS, pH 7.4 with 0.02% Sodium Azide

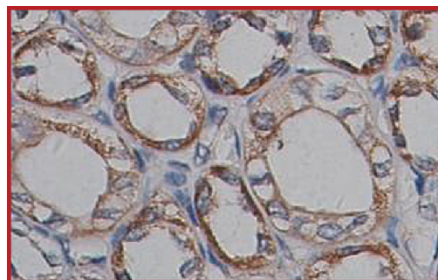
Stability

1yr at -20°C from date of shipment

References

1. Read,R.C., Teare,D.M., Pridmore,A.C., et al. The tumor necrosis factor polymorphism TNF (-308) is associated with susceptibility to meningococcal sepsis, but not with lethality. Crit. Care Med. 37 (4), 1237-1243 (2009).

Sample Data



Immunohistochemical staining of formalin-fixed paraffin-embedded fetal kidney showing membrane staining with Anti-NOTCH 4 at a dilution of 1:100.

Anti-NOTCH 4

Rabbit Polyclonal Antibody

Catalog #	N44-363R
Lot #	Z2014-20
Purification	Affinity chromatography
Stability	1yr at -20°C from date of shipment
Storage & Shipping	Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.