

ASAH2 Polyclonal Antibody

(Catalog # A72338)

Description

ASAH2 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

Q9NR71

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 80-340 of human ASAH2 (NP_001137446.1).

Storage

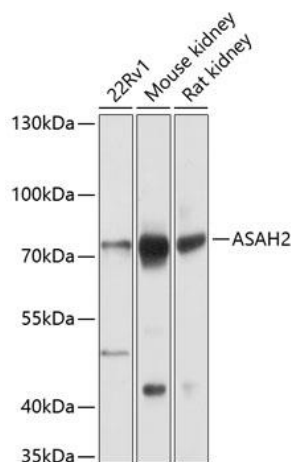
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

ASAH2; BCDase; HNAC1; LCDase; N-CDase; NCDase; neutral ceramidase

Application

WB; Recommended dilution: WB 1:500 - 1:2000



Western blot analysis of extracts of various cell lines, using ASAH2 Polyclonal Antibody at 1:1000 dilution. Secondary Antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25ug per lane. Blocking buffer: 3% nonfat dry milk in TBST. Exposure time: 1s.