



Product Data Sheet

Product Name:	Anti- TMED10 antibody, <i>polyclonal</i> (<i>Transmembrane emp24-like trafficking protein 10</i>)		
Catalog Number:	56015		
Lot Number:	See label on vial		
Product Description:	This antibody is supplied as 50 ug of affinity purified rabbit IgG in 1X PBS (pH 7.4) with 0.05% sodium azide.		
Immunogen:	Rabbit polyclonal anti-TMED10 antibody was raised against a synthetic peptide corresponding to near the C-terminus of human TMP21 protein		
Species Reactivity:	Human, Mouse, Rat		
Validation:	WB:	1: 500-2,000	
	IHC:	1: 400	
	IF:	1: 50	
	Other applications	Not tested	
Storage:	Store at 4 °C for up to one year.		
	For long term storage, aliquot and freeze. Avoid repeated freezing and thawing.		

This product is for *in vitro* research use only.

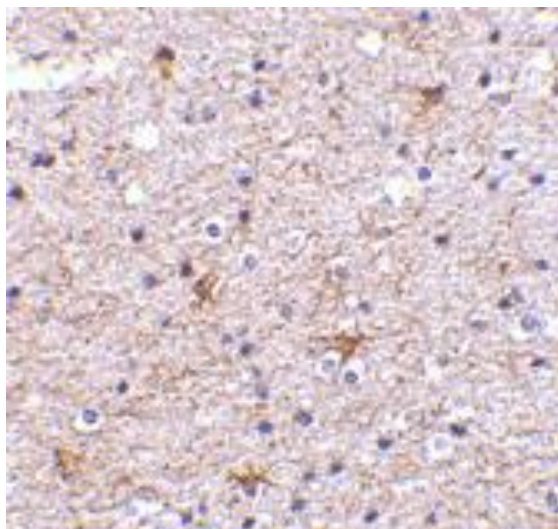
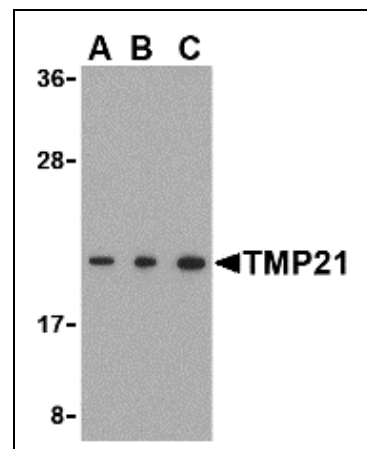


Fig. 1.
IHC of human brain tissue
stained with polyclonal anti-
TMP21 antibody
(Cat#56015)

Fig. 2.
WB of Raji cell lysate
polyclonal anti-TMP21
antibody (Cat# 56015)
(A) 0.5, (B) 1 and (C) 2ug/ml



Target Information:

Official symbol:	TMED10
Official full name:	Transmembrane emp24-like trafficking protein 10
<i>Alternative Names:</i>	21 kDa transmembrane-trafficking protein, S31III125, Tmp-21-I, transmembrane protein Tmp21, p23, p24 family protein delta-1, p24delta
Organism:	Human
Length:	219 amino acids
Predicted MW:	25.0 kDa
Domains:	GOLD domain
Processing:	The sequence is further processed into the mature form.
Subunit structure:	Predominantly homodimeric and to lesser extent monomeric in endoplasmic reticulum.
Tissue expression:	Ubiquitous.
Subcellular location:	Golgi apparatus, cis-Golgi network membrane, trans-Golgi network membrane, single-pass type I membrane protein, melanosome, endoplasmic reticulum membrane, endoplasmic reticulum-Golgi intermediate compartment membrane, cytoplasmic vesicle, secretory vesicle membrane. Note: Identified by mass spectrometry in melanosome fractions from stage I to stage IV. Cycles between compartments of the early secretory pathway.
Function:	Tmp21 is a protein involved in vesicular protein trafficking. Mainly functions in the early secretory pathway. Thought to act as cargo receptor at the luminal side for incorporation of secretory cargo molecules into transport vesicles and to be involved in vesicle coat formation at the cytoplasmic side. In COPII vesicle-mediated anterograde transport involved in the transport of GPI-anchored proteins and proposed to act together with TMED2 as their cargo receptor; the function specifically implies SEC24C and SEC24D of the COPII vesicle coat and lipid raft-like microdomains of the ER. Recognizes GPI anchors structural remodeled in the ER by PGAP1 and MPPE1. In COPI vesicle-mediated retrograde transport involved in the biogenesis of COPI vesicles and vesicle coat recruitment. On Golgi membranes, acts as primary receptor for ARF1-GDP which is involved in COPI-vesicle formation. Increases coatomer-dependent GTPase-activating activity of ARFGAP2. Involved in trafficking of G protein-coupled receptors (GPCRs). Regulates F2LR1, OPRM1 and P2RY4 exocytic trafficking from the Golgi to the plasma membrane thus contributing to receptor resensitization. Involved in trafficking of amyloid beta A4 protein and soluble APP-beta release (independent of modulation of gamma-secretase activity). As part of the presenilin-dependent gamma-

secretase complex regulates gamma-cleavages of the amyloid beta A4 protein to yield amyloid-beta 40 (Abeta40). Involved in organization of the Golgi apparatus.

Disease association: None identified

Source of target information is from HUGO Gene Nomenclature Committee, UniProt, ExPASy PROSITE.