

PRODUCT INFORMATION



Tafazzin Polyclonal Antibody

Item No. 44967

Overview and Properties

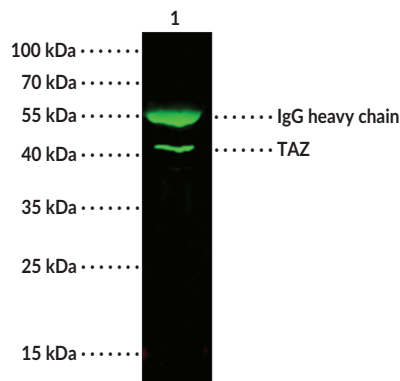
Contents:	This vial contains 100 µl of protein A-affinity purified polyclonal antibody.
Synonyms:	BTHS, CMD3A, EFE2, LVNCX, Protein G4.5, TAZ
Immunogen:	Synthetic peptide corresponding to the C-terminal region of human tafazzin
Cross Reactivity:	(+) Tafazzin
Species Reactivity:	(+) Human
Form:	Liquid
Storage:	-80°C (as supplied)
Stability:	≥1 year
Storage Buffer:	0.2 µm filtered solution in PBS
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry using paraffin-embedded samples (IHC-P), immunoprecipitation (IP), and Western blot (WB); the recommended starting dilution is 1:5,000-1:20,000 for IHC-P, 0.2-1 µl/mg of lysate for IP, and 1:500-1:1,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Lane 1: A549 whole cell lysate (30 µg)

WB of Tafazzin Polyclonal Antibody at a dilution of 1:500.



Lane 1: A549 whole cell lysate (500 µg), Tafazzin Polyclonal Antibody (500 µl), and immunomagnetic beads (60 µg)

WB of Tafazzin Polyclonal Antibody at a dilution of 1:500.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

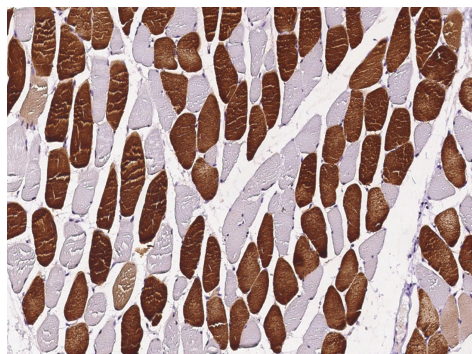
Copyright Cayman Chemical Company, 01/07/2026

CAYMAN CHEMICAL

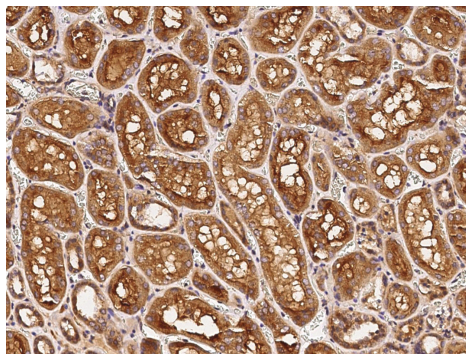
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

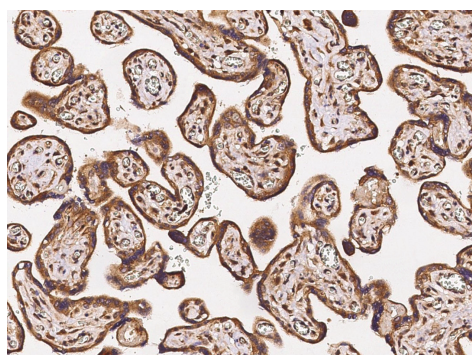
Images cont.



Immunohistochemistry staining of formalin-fixed and paraffin-embedded human TAZ in human skeletal muscle cells using Tafazzin Polyclonal Antibody at a dilution of 1:10,000.



Immunohistochemistry staining of formalin-fixed and paraffin-embedded human TAZ in human kidney cells using Tafazzin Polyclonal Antibody at a dilution of 1:10,000.



Immunohistochemistry staining of formalin-fixed and paraffin-embedded human TAZ in human placenta cells using Tafazzin Polyclonal Antibody at a dilution of 1:10,000.

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Tafazzin is a transacylase involved in cardiolipin remodeling.¹ It is composed of a transmembrane domain, an HX4D motif, two targeting domains flanking the transacylase active site, and a membrane anchor. Tafazzin is expressed in a wide variety of tissues, including cardiac and skeletal muscle, and is localized to the inner and outer mitochondrial membranes.^{1,2} It transfers acyl chains from phospholipids to monolysocardiolipin (MLCL) to form mature cardiolipin, with a preference for transferring linoleic acid from phosphatidylcholine to MLCL.^{1,3} Due to its role in determining cardiolipin composition, tafazzin activity indirectly affects mitochondrial function, including oxidative phosphorylation, oxidative stress repair, and apoptosis.¹ Mutations in *TAFAZZIN* are associated with Barth syndrome, an X-linked disorder characterized by cardiovascular, musculoskeletal, and neurological pathologies.³ Cayman's Tafazzin Polyclonal Antibody can be used for immunohistochemistry using paraffin-embedded samples (IHC-P), immunoprecipitation (IP), and Western blot (WB) applications.

References

1. Garlid, A.O., Schaffer, C.T., Kim, J., *et al.* TAZ encodes tafazzin, a transacylase essential for cardiolipin formation and central to the etiology of Barth syndrome. *Gene* **726**, 144148 (2020).
2. Lu, B., Kelher, M.R., Lee, D.P., *et al.* Complex expression pattern of the Barth syndrome gene product tafazzin in human cell lines and murine tissues. *Biochem. Cell Biol.* **82(5)**, 569-576 (2004).
3. Chin, M.T. and Conway, S.J. Role of Tafazzin in mitochondrial function, development and disease. *J. Dev. Biol.* **8(2)**, 10 (2020).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM