

PRODUCT INFORMATION

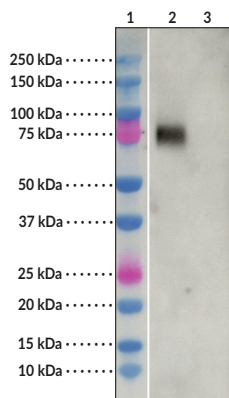
MARCH6 Monoclonal Antibody (Clone 26F12C)

Item No. 44928

Overview and Properties

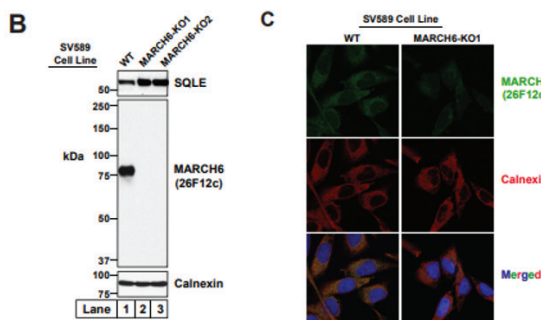
Contents:	This vial contains 100 µg of protein G-purified monoclonal antibody
Synonyms:	Doa10 Homolog, E3 Ubiquitin-protein Ligase MARCH6, MARCHF6, MARCH-VI, Membrane-associated RING-CH Protein VI, Membrane-associated RING-CH-type Finger 6, Membrane-associated RING Finger (C3HC4) 6, Membrane-associated RING Finger Protein 6, RING Finger Protein 176, RNF176, TEB4
Immunogen:	Recombinant human MARCH6 ^{C9A}
Cross Reactivity:	(+) MARCH6
Species Reactivity:	(+) Human; other species not tested
Uniprot No.:	O60337
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
Clone:	26F12C
Host:	Mouse
Isotype:	IgG1K
Applications:	Western Blot (WB); the recommended starting dilution is 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically. Antibody has also been published to be effective for immunofluorescence (IF)

Images



Lane 1: MW Markers
Lane 2: SV-589 Cell Lysate - Wild Type
Lane 3: SV-589 Cell Lysate - Knockout

Western Blot. Reduced samples were loaded into a 4-20% SDS-PAGE gel and transferred to nitrocellulose followed by blocking with 5% NFD (with 0.01% Tween) and probing with MARCH6 Monoclonal Antibody (Clone 26F12C) (Item No. 44928) and Goat Anti-Mouse IgG HRP (Item No. 10004302).



Western blot and immunofluorescence using MARCH6 Monoclonal Antibody (Clone 26F12C) in two CRISPR-generated MARCH6 knockout SV-589 cell lines.

doi: <https://doi.org/10.1016/j.jlr.2024.100650>

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, 06/30/2026

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



Description

Membrane-associated RING-CH-type finger 6 (MARCH6) is an E3 ubiquitin ligase.^{1,2} It is composed of an N-terminal C4HC3-type RING domain, which has catalytic activity, and 14 transmembrane domains. MARCH6 is ubiquitously expressed and localizes to the endoplasmic reticulum (ER).^{1,3} It plays a role in ER-associated degradation (ERAD), where it targets substrates involved in cholesterol synthesis and trafficking, lipid droplet formation, bile salt transport, thyroid hormone activation, and G-protein signaling for proteasomal degradation.^{1,4} MARCH6 is also involved NADPH sensing, ferroptosis, and redox homeostasis.⁵ Intratumoral expression of *MARCHF6* is increased in a variety of cancers, and intronic TTTTA/TTTCA repeat expansions in *MARCHF6* are associated with familial adult myoclonic epilepsy (FAME).^{5,6} Cayman's MARCH6 Monoclonal Antibody (Clone 26F12C) can be used for immunofluorescence (IF) and Western blot (WB) applications. The antibody recognizes MARCH6 at approximately 75 kDa from human samples using 4-20% SDS-PAGE.

References

1. Lin, H., Li, S., and Shu, H.-B. The membrane-associated MARCH E3 ligase family: Emerging roles in immune regulation. *Front. Immunol.* **10**, 1751 (2019).
2. Samji, T., Hong, S., and Means, R.E. The membrane associated RING-CH proteins: A family of E3 ligases with diverse roles through the cell. *Int. Sch. Res. Notices* **2014**, 637295 (2014).
3. Xu, S., Donnelly, L., Kober, D.L., et al. Development of a monoclonal antibody to study MARCH6, an E3 ligase that regulates proteins that control lipid homeostasis. *J. Lipid Res.* **65(11)**, 100650 (2024).
4. Scott, N.A., Sharpe, L.J., and Brown, A.J. The E3 ubiquitin ligase MARCHF6 as a metabolic integrator in cholesterol synthesis and beyond. *Biochim. Biophys. Acta Mol. Cell Biol. Lipids* **1866(1)**, 158837 (2021).
5. Zhen, Y., Chen, S., Niu, B., et al. Unraveling MARCH6's role in cancer progression and metabolism from protein homeostasis to oncogenesis. *Pharmacol. Res.* **229**, 108236 (2026).
6. Florian, R.T., Kraft, F., Leitão, E., et al. Unstable TTTTA/TTTCA expansions in MARCH6 are associated with familial adult myoclonic epilepsy type 3. *Nat. Commun.* **10(1)**, 4919 (2019).

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