

PRODUCT INFORMATION



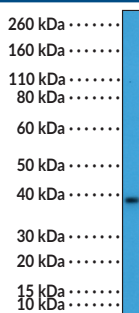
Musashi-2 Rabbit Monoclonal Antibody (Clone RM422)

Item No. 35315

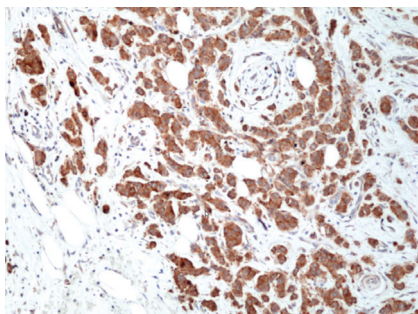
Overview and Properties

Contents:	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
Synonyms:	MSI2, RNA-Binding Protein Musashi Homolog 2
Immunogen:	Peptide from the internal region of Musashi-2
Cross Reactivity:	(+) Musashi-2
Species Reactivity:	(+) Human, mouse
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS, with 50% glycerol, 1% BSA, and 0.09% sodium azide
Clone:	RM422
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:100-1:400 for IHC and 1:100-1:2,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

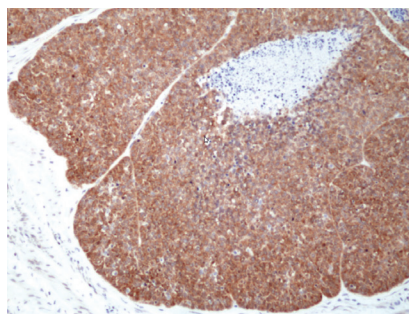
Images



WB of mouse brain tissue lysate using Musashi-2 Rabbit Monoclonal Antibody (Clone RM422) at a dilution of 1:2,000.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human breast ductal carcinoma tissue using Musashi-2 Rabbit Monoclonal Antibody (Clone RM422) at a dilution of 1:400.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human esophageal squamous carcinoma tissue using Musashi-2 Rabbit Monoclonal Antibody (Clone RM422) at a dilution of 1:100.

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, 01/19/2024

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

Musashi-2 is an RNA-binding protein that regulates target mRNA stability and translation.^{1,2} It contains two RNA recognition motifs at its N-terminus that each contain two highly conserved core motifs, RNP-1 and RNP-2, and a C-terminal domain that is subject to alternative splicing to produce a short or long isoform of Musashi-2.^{3,4} Musashi-2 is ubiquitously expressed and localizes to the nucleus and cytoplasm.³⁻⁵ It is involved in the differentiation and self-renewal of stem cells, including embryonic, neural, and hematopoietic stem cells, as well as in the migration and proliferation of keratinocytes.^{1,2} Knockdown of *MSI2*, the gene encoding Musashi-2, induces cell cycle arrest at the G₁ phase and apoptosis in keratinocytes.² Chromosomal translocations leading to the fusion of *MSI2* with *HOXA9*, resulting in expression of Musashi-2-HOXA9 fusion proteins, are associated with chronic myeloid leukemia (CML).⁴ Cayman's Musashi-2 Rabbit Monoclonal Antibody (Clone RM422) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

References

1. Sutherland, J.M., McLaughlin, E.A., Hime, G.R., *et al.* The Musashi family of RNA binding proteins: Master regulators of multiple stem cell populations. *Transcriptional and translational regulation of stem cells*. Hime, G. and Abud, H., editors, *Springer* (2013).
2. Bennett, C.G., Riemondy, K., Chapnick, D.A., *et al.* Genome-wide analysis of Musashi-2 targets reveals novel functions in governing epithelial cell migration. *Nucleic Acids Res.* **44(8)**, 3788-3800 (2016).
3. Sakakibara, S., Nakamura, Y., Satoh, H., *et al.* RNA-binding protein Musashi2: Developmentally regulated expression in neural precursor cells and subpopulations of neurons in mammalian CNS. *J. Neurosci.* **21(20)**, 8091-8107 (2001).
4. Barbouti, A., Höglund, M., Johansson, B., *et al.* A novel gene, *MSI2*, encoding a putative RNA-binding protein is recurrently rearranged at disease progression of chronic myeloid leukemia and forms a fusion gene with *HOXA9* as a result of the cryptic t(7;17)(p15;q23). *Cancer Res.* **63(6)**, 1202-1206 (2003).
5. Sutherland, J.M., Sobinoff, A.P., Gunter, K.M., *et al.* Knockout of RNA binding protein *MSI2* impairs follicle development in the mouse ovary: Characterization of *MSI1* and *MSI2* during folliculogenesis. *Biomolecules* **5(3)**, 1228-1244 (2015).

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