

# PRODUCT INFORMATION



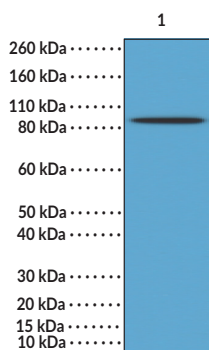
## SATB2 (C-Term) Rabbit Monoclonal Antibody (Clone RM365)

Item No. 32290

### Overview and Properties

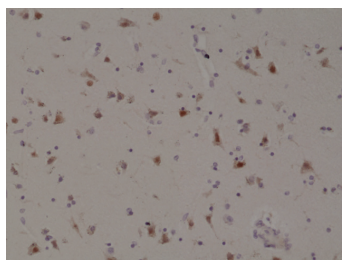
<b>Contents:</b>	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
<b>Synonyms:</b>	Special AT-rich Sequence-Binding Protein 2
<b>Immunogen:</b>	Peptide from the C-terminal region of human DNA-binding protein SATB2
<b>Cross Reactivity:</b>	(+) SATB2
<b>Species Reactivity:</b>	(+) Human
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Storage Buffer:</b>	PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide
<b>Clone:</b>	RM365
<b>Host:</b>	Rabbit
<b>Isotype:</b>	IgG
<b>Applications:</b>	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:200-1:1,000 for IHC and 1:1,000-1:2,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Images

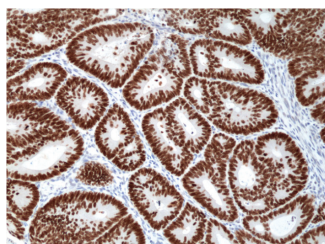


Lane 1: K562 cell lysate

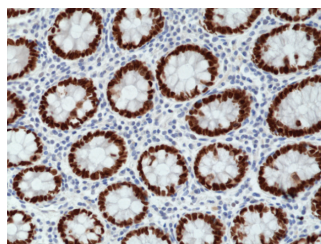
WB of K562 cell lysate using SATB2 (C-Term) Rabbit Monoclonal Antibody (Clone RM365) at dilution of 1:1,000.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human brain tissue using SATB2 (C-Term) Rabbit Monoclonal Antibody (Clone RM365) at a dilution of 1:200.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human colon cancer tissue using SATB2 (C-Term) Rabbit Monoclonal Antibody (Clone RM365) at a dilution of 1:1,000.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human colon tissue using SATB2 (C-Term) Rabbit Monoclonal Antibody (Clone RM365) at a dilution of 1:1,000.

**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, 02/15/2024

**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

---

Special AT-rich sequence-binding protein 2 (SATB2) is a transcription factor.<sup>1</sup> It is composed of two CUT domains and a homeobox domain, which bind DNA matrix attachment regions (MARs) and promote chromatin remodeling.<sup>1,2</sup> SATB2 is expressed in superficial cortical layers during embryonic development, where it regulates the formation of callosal projection neurons by repressing *Ctip2* expression, and by hippocampal and cortical pyramidal neurons in the adult brain, where it has roles in synaptic plasticity and neurotransmission.<sup>2,3</sup> It has additional functions in B cell and osteoblast differentiation.<sup>2</sup> SATB2 has been used as a marker for primary or metastatic colorectal cancer.<sup>4</sup> SATB2 variants have been found in patients with SATB2-associated syndrome, an autosomal dominant disorder characterized by intellectual disability and craniofacial abnormalities, as well as in patients with schizophrenia.<sup>1,5</sup> Cayman's SATB2 (C-Term) Rabbit Monoclonal Antibody (Clone RM365) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

## References

---

1. Zarate, Y.A., Bosanko, K.A., Caffrey, A.R., *et al.* Mutation update for the SATB2 gene. *Hum. Mutat.* **40(8)**, 1013-1029 (2019).
2. Alcamo, E.A., Chirivella, L., Dautzenberg, M., *et al.* Satb2 regulates callosal projection neuron identity in the developing cerebral cortex. *Neuron* **57(3)**, 364-377 (2008).
3. Cera, I., Whitton, L., Donohoe, G., *et al.* Genes encoding SATB2-interacting proteins in adult cerebral cortex contribute to human cognitive ability. *PLoS Genet.* **15(2)**, e1007890 (2019).
4. Zhang, Y.-J., Chen, J.-W., He, X.-S., *et al.* SATB2 is a promising biomarker for identifying a colorectal origin for liver metastatic adenocarcinomas. *EBioMedicine* **28**, 62-69 (2018).
5. Zhou, J., Chen, J., Xu, W., *et al.* Common variants in SATB2 are associated with schizophrenia in Uyghur Chinese population. *Psychiatr. Genet.* **29(4)**, 120-126 (2018).

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