

PRODUCT INFORMATION



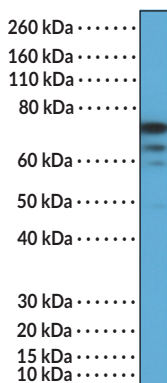
FOXP1 (C-Term) Rabbit Monoclonal Antibody (RM402)

Item No. 32328

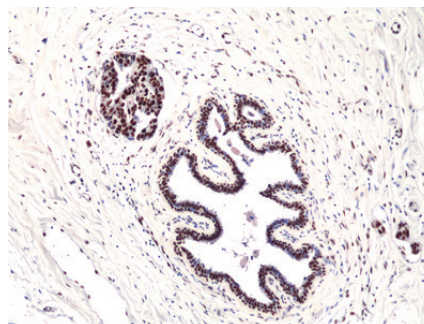
Overview and Properties

Contents:	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
Synonyms:	Forkhead Box P1, Glutamine-rich Factor 1, QRF1
Immunogen:	Peptide from the C-terminal region of human FOXP1
Cross Reactivity:	(+) FOXP1
Species Reactivity:	(+) Human
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide
Clone:	RM402
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:100-1:200 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



WB of 293 cell lysate using FOXP1 (C-Term) Rabbit Monoclonal Antibody (RM402) at a dilution of 1:1,000.



Immunohistochemical staining of formalin-fixed and paraffin-embedded breast ductal carcinoma tissue using FOXP1 (C-Term) Rabbit Monoclonal Antibody (RM402) at a 1:200 dilution.

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, 02/21/2024

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

Forkhead box P1 (FOXP1) is a member of the FOX family of transcription factors and a transcriptional repressor with roles in cardiac valve morphogenesis, B and T cell development, monocyte differentiation, and lung development.¹ It is composed of an N-terminal polyglutamine tract, a zinc finger domain, a leucine zipper, and a C-terminal forkhead domain.^{1,2} FoxP1 is ubiquitously expressed and localized to the nucleus where it binds to DNA to regulate transcription. Increased expression of *FOXP1* is associated with poor prognosis in patients with diffuse large B cell lymphoma (DLBCL).³ Mutations in *FOXP1* are associated with language impairment, speech delay, and grammar defects in patients with autism spectrum disorder.^{2,4} Cayman's FOXP1 (C-Term) Rabbit Monoclonal Antibody (RM402) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

References

1. Koon, H.B., Ippolito, G.C., Banham, A.H., *et al.* FOXP1: A potential therapeutic target in cancer. *Expert Opin. Ther. Targets* **11(7)**, 955-965 (2007).
2. Bowers, J.M., and Konopka, G. The role of the FOXP family of transcription factors in ASD. *Dis. Markers* **33(5)**, 251-260 (2012).
3. Tzankov, A., Zlobec, I., Went, P., *et al.* Prognostic immunophenotypic biomarker studies in diffuse large B cell lymphoma with special emphasis on rational determination of cut-off scores. *Leuk. Lymphoma*. **51(2)**, 199-212 (2010).
4. Bacon, C. and Rappold, G.A. The distinct and overlapping phenotypic spectra of FOXP1 and FOXP2 in cognitive disorders. *Hum. Genet.* **131(11)**, 1687-1698 (2012).

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