

PRODUCT INFORMATION

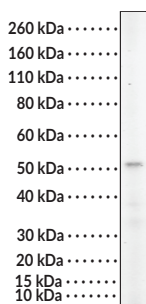
GATA3 Rabbit Monoclonal Antibody (Clone RM482)

Item No. 42776

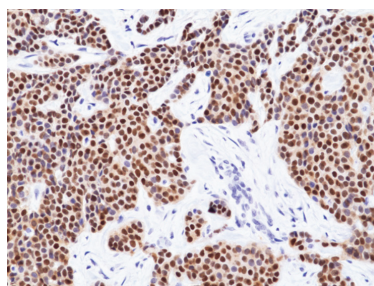
Overview and Properties

Contents:	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
Synonyms:	GATA-binding Factor 3, Trans-acting T cell-specific Transcription Factor GATA3
Immunogen:	Peptide corresponding to human GATA3
Cross Reactivity:	(+) GATA3
Species Reactivity:	(+) Human, mouse, rat
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide
Clone:	RM482
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:100-1:200 for IHC and 1:100-1:500 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

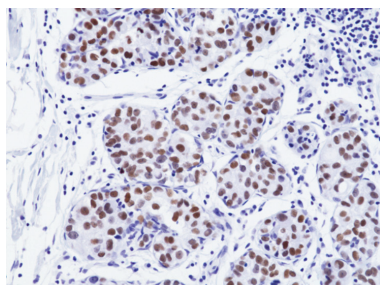
Images



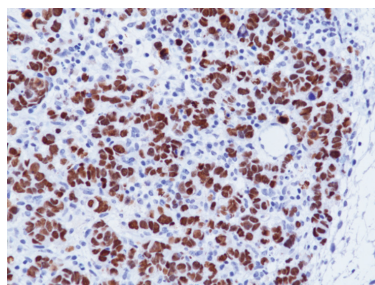
WB of Jurkat cell lysates using GATA3 Rabbit Monoclonal Antibody (Clone RM482) at a dilution of 1:500.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human breast ductal carcinoma tissue using GATA3 Rabbit Monoclonal Antibody (Clone RM482) at a dilution of 1:100 for one hour at room temperature.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human breast pleomorphic lobular carcinoma tissue using GATA3 Rabbit Monoclonal Antibody (Clone RM482) at a dilution of 1:100 for one hour at room temperature.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human metastatic breast cancer in ovary tissue using GATA3 Rabbit Monoclonal Antibody (Clone RM482) at a dilution of 1:100 for one hour at room temperature.

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, 03/17/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

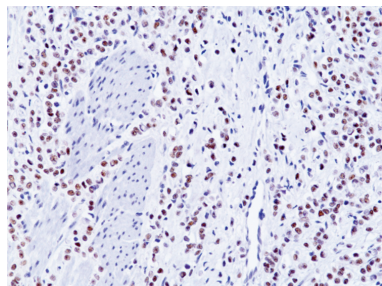
PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

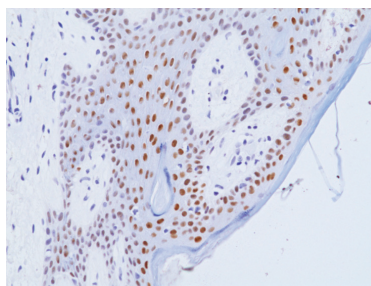
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

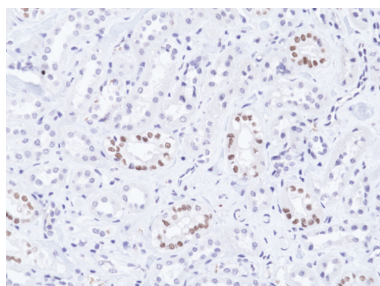
Images Cont.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human metastatic breast cancer in stomach tissue using GATA3 Rabbit Monoclonal Antibody (Clone RM482) at a dilution of 1:100 for one hour at room temperature.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human normal skin tissue using GATA3 Rabbit Monoclonal Antibody (Clone RM482) at a dilution of 1:100 for one hour at room temperature.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human normal kidney tissue using GATA3 Rabbit Monoclonal Antibody (Clone RM482) at a dilution of 1:100 for one hour at room temperature.

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

GATA3 is a zinc finger protein and transcription factor with roles in cell differentiation and organogenesis.^{1,2} It is composed of two N-terminal transactivating domains and two C-terminal zinc finger domains.³ GATA3 is expressed in hematopoietic cells, neurons, adipose tissue, mammary glands, renal tissue, skin, hair, parathyroid glands, and skeletal muscle.¹ It functions as a primary transcription factor in germ cell specification programs and regulates the redistribution of T cell populations from Th2 to Th1 T helper cells. Mutations in GATA3 are associated with the autosomal dominant disorder hypoparathyroidism, sensorineural deafness, and renal dysplasia (HDR) syndrome, also known as Barakat syndrome.³ GATA3 downregulation is associated with poor prognosis and increased mortality in patients with urothelial malignancies.² Cayman's GATA3 Rabbit Monoclonal Antibody (Clone RM482) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

References

1. Bacha, R., Pedersen, S., Ismail, R., *et al.* GATA3: Orchestrating cellular fate through differentiation and proliferation. *Biochim. Biophys. Acta Mol. Cell Res.* **1872**(2), 119893 (2025).
2. Khazaeli Najafabadi, M., Mirzaeian, E., Memar Montazerin, S., *et al.* Role of GATA3 in tumor diagnosis: A review. *Pathol. Res. Pract.* **226**:153611, (2021).
3. Huang, B., Li, S., Chai, Y., *et al.* A novel GATA3 frameshift mutation causes hypoparathyroidism, sensorineural deafness, and renal dysplasia syndrome. *Mol. Genet. Metab. Rep.* **38**:101063, (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